

Quantikine™ IVDELISA

Human Erythropoietin

Quantikine™ IVD ELISA

Human Erythropoietin Immunoassay

REF DEP00B

This package insert must be read in its entirety before using this product.

IVD For *In Vitro* Diagnostic use.

TABLE OF CONTENTS

SECTION	PAGE
NAME AND INTENDED USE.....	1
SUMMARY AND EXPLANATION.....	1
ASSAY PRINCIPLE.....	2
LIMITATIONS.....	2
WARNINGS AND PRECAUTIONS.....	3
INDICATIONS OF INSTABILITY OR DETERIORATION.....	3
OTHER SUPPLIES REQUIRED.....	4
INSTRUMENTS.....	4
MATERIALS PROVIDED & STORAGE CONDITIONS.....	5
SPECIMEN COLLECTION AND STORAGE.....	6
REAGENT PREPARATION.....	6
DILUTION OF SPECIMENS WITH HIGH EPO CONCENTRATIONS.....	6
ASSAY PROCEDURE.....	7
ASSAY PROCEDURE SUMMARY.....	8
CALCULATION OF RESULTS.....	9
TYPICAL DATA.....	9
QUALITY CONTROL.....	10
CALIBRATION.....	10
EXPECTED VALUES.....	11
PERFORMANCE CHARACTERISTICS.....	11
TROUBLESHOOTING GUIDE.....	15
REFERENCES.....	16
PLATE LAYOUT.....	17

 **MANUFACTURED AND DISTRIBUTED BY:**

USA R&D Systems, Inc.

614 McKinley Place NE, Minneapolis, MN 55413

TEL: 800 343 7475 612 379 2956

FAX: 612 656 4400

E-MAIL: info@bio-techne.com

NAME AND INTENDED USE

Quantikine™ IVD Human Erythropoietin ELISA **R&D Systems™ Inc. Catalog Number DEP00B**

Enzyme linked immunosorbent assay (ELISA) for the quantitative determination of Erythropoietin (Epo) concentration in human serum and plasma as an aid in the diagnosis of anemia and polycythemia.

SUMMARY AND EXPLANATION

Erythropoietin, a glycoprotein (~30,400 Daltons) produced primarily by the kidney, is the principal factor regulating red blood cell production (erythropoiesis) in mammals. Renal production of Epo is regulated by changes in oxygen availability. Under conditions of hypoxia, the level of Epo in the circulation increases and this leads to increased production of red blood cells.

The over-expression of Epo may be associated with certain pathophysiological conditions (1, 2). Polycythemia exists when there is an overproduction of red blood cells (RBCs). Primary polycythemia, such as polycythemia vera, are caused by Epo-independent growth of erythrocytic progenitors from abnormal stem cells and low to normal levels of Epo are found in the serum of affected patients. On the other hand, various types of secondary polycythemia are associated with the production of higher than normal levels of Epo. The overproduction of Epo may be an adaptive response associated with conditions that produce tissue hypoxia, such as living at high altitude, chronic obstructive pulmonary disease, cyanotic heart disease, sleep apnea, high-affinity hemoglobinopathy, smoking, or localized renal hypoxia (1, 2). In other instances, excessive Epo levels are the result of production by neoplastic cells. Cases of increased Epo production and erythrocytosis have been reported for patients with renal carcinomas (3), benign renal tumors (4), Wilms' tumors, hepatomas (5), liver carcinomas (6), cerebellar hemangioblastomas (3, 7, 8), adrenal gland tumors (9), smooth muscle tumors (3, 9), and leiomyomas (10).

Deficient Epo production is found in conjunction with certain forms of anemias. These include anemia of renal failure and end-stage renal disease (1, 2, 11), anemias of chronic disorders [chronic infections (1), autoimmune diseases (1), rheumatoid arthritis (12), AIDS (13), malignancies (14)], anemia of prematurity (2), anemia of hypothyroidism (2), and anemia of malnutrition (2). Many of these conditions are associated with the generation of IL-1 and TNF- α , factors that have been shown to be inhibitors of Epo activity (1, 15). Other forms of anemias, on the other hand, are due to Epo-independent causes and affected individuals show elevated levels of Epo (2). These forms include aplastic anemias, iron deficiency anemias, thalassemias, megaloblastic anemias, pure red cell aplasias, and myelodysplastic syndromes.

Methods for determining Epo concentration in serum have been based historically on the ^{59}Fe exhypoxic polycythemic mouse *in vivo* bioassay (16). The Quantikine™ IVD Human Epo ELISA uses a monoclonal antibody and polyclonal antibody conjugate in a sandwich ELISA format to provide a method that is quicker, more sensitive and more specific than the bioassay. The assay is designed to measure Epo levels in serum or plasma in less than 4.5 hours, or less than 2.5 hours using the shaker protocol.

ASSAY PRINCIPLE

The Quantikine™ IVD Human Epo ELISA is based on the double-antibody sandwich method. Microplate wells, precoated with a monoclonal antibody specific for Epo are incubated with specimen or standard. Erythropoietin binds to the immobilized antibody on the plate. After removing excess specimen or standard, wells are incubated with an anti-Epo polyclonal antibody conjugated to horseradish peroxidase. During the second incubation, the antibody-enzyme conjugate binds to the immobilized Epo. Excess conjugate is removed by washing. A chromogen is added to the wells and is oxidized by the enzyme reaction to form a blue colored complex. The reaction is stopped by the addition of acid, which turns the blue to yellow. The amount of color generated is directly proportional to the amount of conjugate bound to the Epo antibody complex, which, in turn, is directly proportional to the amount of Epo in the specimen or standard. The absorbance of this complex is measured and a standard curve is generated by plotting absorbance versus the concentration of the Epo standards. The Epo concentration of the unknown specimen is determined by comparing the optical density of the specimen to the standard curve. The standards used in this assay are recombinant human Epo calibrated against the Third International Standard (11/170), a recombinant form of human Erythropoietin.

LIMITATIONS

- The results of this assay should be used in conjunction with information available from clinical evaluations and other diagnostic procedures.
- No drugs have been investigated for assay interference.
- If specimens generate values higher than the highest standard, dilute the specimens in Specimen Diluent and repeat the assay.
- Any variation in operator, pipetting technique, washing technique, incubation time or temperature, and kit age can cause variation in binding.

WARNINGS AND PRECAUTIONS

IVD For *In Vitro* Diagnostic Use

- Do not use kit reagents beyond the kit expiration date.
 - For best results, each laboratory should validate a specific assay method (benchtop or shaker) and perform all assays by that method.
 - In order to minimize within assay variation, it is recommended that the assay be pipetted within 20 minutes.
 - Do not substitute kit reagents with those from other lots or other sources.
 - Do not expose kit reagents to strong light during storage or incubation.
 - Avoid contact of kit reagents with oxidizing agents and metal.
 - Exposure to sodium azide will inactivate the conjugate.
 - Do not pipette by mouth.
 - Do not smoke or eat in areas where kit reagents or specimens are handled.
 - Avoid contact of skin and mucous membranes with kit reagents or specimens.
 - If any reagents come in contact with eyes, skin, or mucous membranes, wash with copious amounts of water and contact a physician.
 - Handle all serum and materials in contact with serum in accordance with CLSI guidelines for preventing the transmission of blood-borne pathogens during laboratory procedures.
 - Incubation times and temperatures other than those specified may yield erroneous results.
 - Contamination of kit reagents may yield erroneous results.
 - If possible, use plastic disposable pipettes, tips and containers for reagent preparation and storage. Glassware used should be rinsed thoroughly with 1 N sulfuric or 1 N hydrochloric acid followed by at least three washes of deionized water. No acid or detergent residues should remain on the glassware.
 - Use polypropylene or high density polyethylene (HDPE) test tubes for specimen dilutions, if needed. **DO NOT USE GLASS TUBES.**
 - Erythropoietin Assay Diluent contains sodium azide which may react with lead and copper plumbing to form explosive metal azides. Flush with large volumes of water during disposal.
- Note:** Users shall report any serious incident that has occurred in relation to the device to R&D Systems™, Inc. at 1-800-343-7475 (USA and Canada) or 1-612-379-2956 and the competent authority of the Member State in which the user and/or patient is established.

INDICATIONS OF INSTABILITY OR DETERIORATION

Precipitates in the reagent solutions are generally considered indications of instability or deterioration. If any of these indications of instability or deterioration are observed, or if the correlation coefficient of the standard curve is less than 0.95, retain the reagent(s) in question at 2-8 °C and contact R&D Systems™, Inc. at 1-800-343-7475 (USA and Canada) or 1-612-379-2956.

OTHER SUPPLIES REQUIRED

- Pipettes and pipette tips
- Horizontal orbital microplate shaker (0.12" orbit) capable of maintaining a speed of 500 ± 50 rpm (required for the Shaker Protocol)
- 20 mL and 500 mL graduated cylinders
- Squirt bottle, manifold dispenser, or automated microplate washer
- Absorbent pad or paper towels for blotting the wells
- Microplate reader capable of measuring absorbance at 450 nm, with the correction wavelength set at 600 nm
- Immunoassay reagent trays
- Deionized or distilled water
- A computer capable of 4 parameter logistic curve fitting for data reduction
- **Erythropoietin Serum Controls** - e.g., Quantikine™ IVD Human Serum Control 1 & 2 and Control 3 (R&D Systems™, Catalog # CEP01: 5 vials each of Levels 1 & 2; Catalog # CEP03: 10 vials of Level 3), or equivalent

INSTRUMENTS

Assay results are measured spectrophotometrically at 450 nm using a microplate reader. For best results, a reference wavelength of 600 nm (540 nm, 570 nm, and 650 nm may also be used) should be included in the measurement to correct for optical imperfections in the polystyrene microplate. Instruments without reference filters may be used, but assay precision may decrease. A microplate reader with an optical density range of 0-3 Optical Density (O.D.) and an accuracy of ± 0.005 O.D. is recommended. Microplate readers with an O.D. range less than 0-3 O.D. may be used, but the range of the assay may be reduced.

MATERIALS PROVIDED & STORAGE CONDITIONS

Store the unopened kit at 2-8 °C. Do not use past kit expiration date.

			DESCRIPTION	STORAGE OF OPENED/DILUTED MATERIAL
SORB			Epo 96 Well Microplate (Part 891647) - 96 well polystyrene microplate (12 strips of 8 wells) coated with a monoclonal antibody specific for recombinant human Epo.	Return unused wells to the foil pouch containing the desiccant pack. Reseal along entire edge of the zip-seal. May be stored for up to 1 month at 2-8 °C.*
CONJ			Epo Conjugate (Part 899576) - 21.5 mL of a polyclonal antibody specific for recombinant human Epo, conjugated to horseradish peroxidase with preservatives.	May be stored for up to 1 month at 2-8 °C.*
CAL	0		Epo 0 mIU/mL Standard (Part 892950) - 2.1 mL of a buffered protein base with preservatives.	
CAL	2.5		Epo 2.5 mIU/mL Standard (Part 893919) - 2.1 mL of recombinant human Epo in a buffered protein base with preservatives.	
CAL	5		Epo 5 mIU/mL Standard (Part 895228) - 2.1 mL of recombinant human Epo in a buffered protein base with preservatives.	
CAL	20		Epo 20 mIU/mL Standard (Part 895230) - 2.1 mL of recombinant human Epo in a buffered protein base with preservatives.	
CAL	50		Epo 50 mIU/mL Standard (Part 895233) - 2.1 mL of recombinant human Epo in a buffered protein base with preservatives.	
CAL	100		Epo 100 mIU/mL Standard (Part 895243) - 2.1 mL of recombinant human Epo in a buffered protein base with preservatives.	
CAL	200		Epo 200 mIU/mL Standard (Part 895259) - 2.1 mL of recombinant human Epo in a buffered protein base with preservatives.	
DIL	AS		Epo Assay Diluent (Part 895217) - 11 mL of a buffered protein base with preservatives. Contains sodium azide.	
DIL	SPE		Epo Specimen Diluent (Part 895218) - 26 mL of a protein stabilized buffer with preservatives.	
WASH	BUF	25X	Epo Wash Buffer Concentrate (Part 895221) - 21 mL of a 25-fold concentrate with preservative.	Room Temperature
TMB	SUBS		Epo TMB Substrate (Part 642881) - 17 mL of tetramethylbenzidine.	
STOP	SOLN		Epo Stop Solution (Part 642879) - 17 mL of a methanesulfonic acid. Caution: <i>The Stop Solution with this kit is an acid solution. Wear protective gloves, clothing, eye, and face protection. Wash hands thoroughly after handling.</i>	
			Plate Covers - 4 adhesive strips.	

* Provided this is within the expiration date of the kit.

SPECIMEN COLLECTION AND STORAGE

Serum - Use a serum separator tube (SST) and allow samples to clot for 30 minutes at room temperature (20-25 °C) before centrifugation for 15 minutes at 1000 x g*. Remove serum and assay immediately or aliquot and store samples at ≤ -20 °C. **Avoid repeated freeze-thaw cycles.**

Plasma - Collect plasma using EDTA as an anticoagulant. Centrifuge for 15 minutes at 1000 x g* within 30 minutes of collection. Assay immediately or aliquot and store samples at ≤ -20 °C. **Avoid repeated freeze-thaw cycles.**

It is recommended that each laboratory standardize the assay using either serum or EDTA plasma specimens.

Lipemic, grossly hemolyzed, or contaminated specimens may yield inaccurate results and should not be tested with this procedure. Further, no drugs have been investigated for assay interference.

Refer to CLSI Guideline: *Procedures for the Handling and Processing of Blood Specimens for Common Laboratory Tests*, (CLSI Document GP44; 4th Edition)

* $g = (1.118 \times 10^{-5}) (\text{radius in cm}) (\text{rpm})^2$

REAGENT PREPARATION

Bring all reagents to room temperature (20-25 °C) before use.

Wash Buffer (1X) - If crystals have formed in the concentrate, warm to room temperature and mix gently until the crystals have completely dissolved. Add 20 mL of Wash Buffer Concentrate to 480 mL of deionized or distilled water to prepare 500 mL of Wash Buffer.

DILUTION OF SPECIMENS WITH HIGH EPO CONCENTRATIONS

If a serum or plasma specimen is above 200 mIU/mL, dilute it with the Epo Specimen Diluent. For example:

- For specimens with Epo concentrations between 200 mIU/mL and 2000 mIU/mL, a 10-fold dilution of the specimen is necessary. Dilute 25 µL of specimen with 225 µL of the Epo Specimen Diluent.
- For specimens with Epo concentrations in excess of 2000 mIU/mL, a higher dilution will be necessary to bring them within the range of the standard curve (*i.e.* 20-fold, 40-fold, *etc.* dilution).

Note: Use polypropylene or high density polyethylene (HDPE) test tubes for specimen dilutions. **DO NOT USE GLASS TUBES.** The use of glass tubes will cause erroneous results due to adsorption of Epo to the glass.

To determine the Epo concentration of the serum or plasma specimen, multiply the result obtained by the dilution factor.

ASSAY PROCEDURE

Bring all reagents and specimens to room temperature (20-25 °C) before use. It is recommended that all standards, controls, and specimens be assayed in duplicate. Benchtop and shaker protocols are provided. The same protocol must be used throughout the assay.

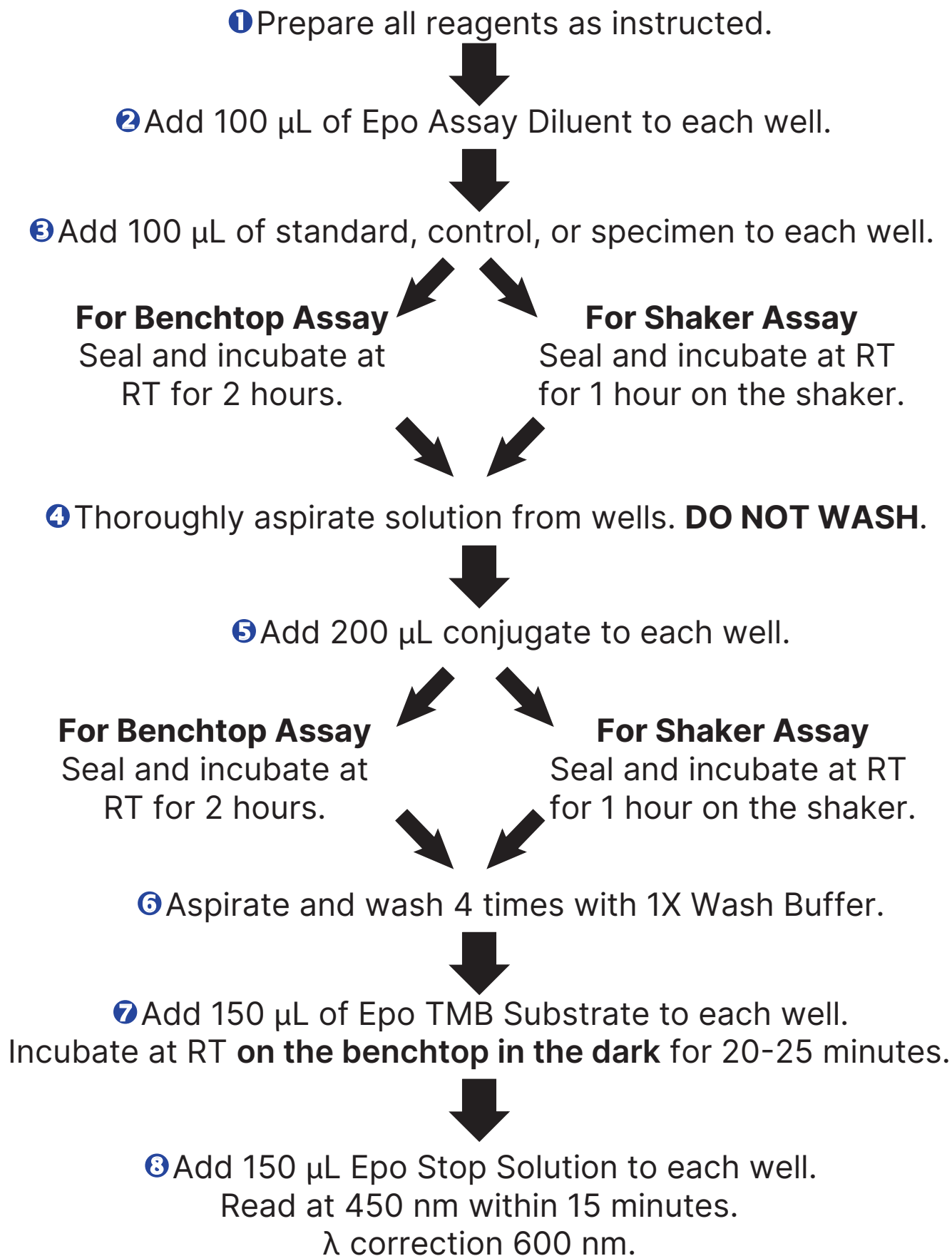
1. Prepare all reagents as directed in the previous section.
2. Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, and reseal.
3. Pipette 100 µL of Epo Assay Diluent into each well.
4. Add 100 µL of standard, control, or specimen per well. Gently tap the plate frame for approximately 1 minute to mix the well contents. Cover the plate with the adhesive strip provided. A plate layout containing a sample diagram of standards, controls, and specimens is shown on page 17.

For Benchtop Protocol: Incubate for 2 hours $\pm$ 5 minutes at room temperature.

For Shaker Protocol: Incubate for 1 hour $\pm$ 5 minutes at room temperature on a horizontal orbital microplate shaker (0.12" orbit) set at 500 $\pm$ 50 rpm.

5. Thoroughly aspirate or decant the contents from each well. Blot dry on clean paper toweling. **Do Not Wash.**
6. Add 200 µL of Epo Conjugate to each well. Cover the plate with a new adhesive strip.
For Benchtop Protocol: Incubate for 2 hours $\pm$ 5 minutes at room temperature.
For Shaker Protocol: Incubate for 1 hour $\pm$ 5 minutes at room temperature on a horizontal orbital microplate shaker.
7. Aspirate each well and wash, repeating the process three times for a total of 4 washes. Wash by filling each well with Wash Buffer (300-400 µL) using a squirt bottle, manifold dispenser, or autowasher. Complete removal of liquid at each step is essential for good performance. After the last wash, remove any remaining Wash Buffer by aspirating or decanting. Invert the plate and blot it against clean paper towels.
8. Add 150 µL of Epo TMB Substrate to each well. Incubate for 20-25 minutes at room temperature **on the benchtop. Protect from light.**
9. Add 150 µL of Epo Stop Solution to each well. If color change does not appear uniform, gently tap the plate to ensure thorough mixing.
10. Determine the optical density (O.D.) of each well within 15 minutes, using a microplate reader set to 450 nm. If wavelength correction is available, set to 600 nm. If wavelength correction is not available, subtract readings at 600 nm from the readings at 450 nm. This subtraction will correct for optical imperfections in the plate. Readings made directly at 450 nm without correction may be higher and less accurate.

ASSAY PROCEDURE SUMMARY



CALCULATION OF RESULTS

Read the absorbance of each well on a microplate reader using 450 nm as the primary wavelength and 600 nm as the reference wavelength (540, 570, or 650 nm is acceptable).

Average the duplicate readings for each standard, control, and specimen and subtract the average 0 mIU/mL standard optical density.

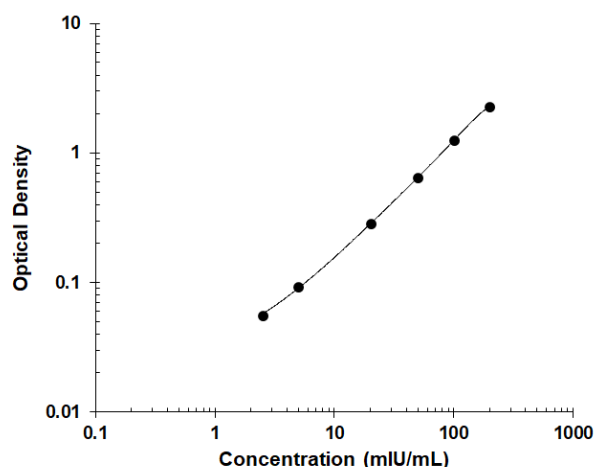
Create a standard curve by reducing the data using software capable of generating a four parameter logistic (4PL) curve-fit. As an alternative, construct a standard curve by plotting the mean absorbance for each standard on the y axis against the concentration on the x axis and draw a best fit curve through the points on the graph. The data may be linearized by plotting the log of the Epo concentrations versus the log of the O.D., and the best fit line can be determined by regression analysis. This procedure will produce an adequate but less precise fit of the data.

Report values for each unknown that reads within the range (2.5-200 mIU/mL) of the assay. For unknown values above the range, see Dilution of Specimens with High Epo Concentrations section. For values below the range, report as undetectable or < 2.5 mIU/mL.

TYPICAL DATA

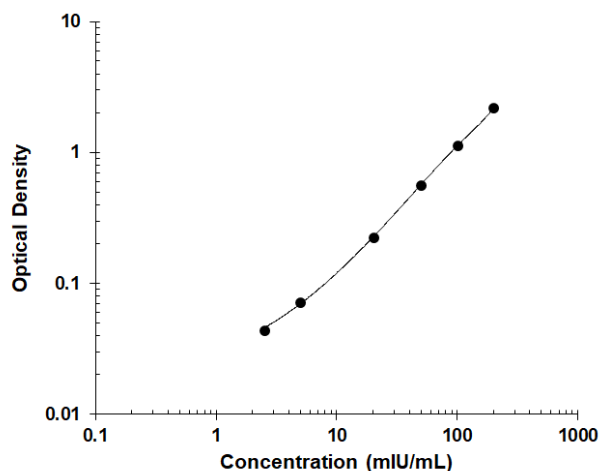
These standard curves are provided for demonstration only. A standard curve should be generated for each set of specimens assayed.

BENCHTOP PROTOCOL



(mIU/mL)	O.D.	Average	Corrected
0	0.008	0.008	
2.5	0.063	0.064	0.056
5.0	0.098	0.100	0.092
20	0.289	0.291	0.283
50	0.655	0.656	0.648
100	1.247	1.258	1.250
200	2.278	2.288	2.280

SHAKER PROTOCOL



(mIU/mL)	O.D.	Average	Corrected
0	0.010	0.012	—
2.5	0.054	0.055	0.044
5.0	0.083	0.084	0.072
20	0.235	0.236	0.225
50	0.575	0.579	0.568
100	1.121	1.137	1.125
200	2.226	2.226	2.215

QUALITY CONTROL

Each testing laboratory should establish a quality control program to monitor the performance of the Quantikine™ IVD Human Epo Immunoassay. As part of this program, controls with known Epo concentrations (available from R&D Systems™) should be run in each assay. R&D Systems™ recommends that at least two controls be run to verify the performance of the assay method. A control in the mid to upper end of the normal range and a control in the mid-region of the assay curve are good choices for the day to day evaluation of the performance of the assay. A control may also be placed at the upper end of the assay curve to evaluate the performance of the upper end of the assay. If the values obtained are not within their established ranges, the assay results may be invalid.

The results of an individual assay are valid if the control values obtained are within the published ranges of a commercially available control or the established range for an in-house control. The correlation coefficient of the fitted standard curve should be ≥ 0.95 .

CALIBRATION

This immunoassay is calibrated against a highly purified recombinant human Epo protein produced at R&D Systems™.

The NIBSC/WHO Human Erythropoietin International Standard 11/170 (recombinant human EPO) was evaluated in this kit. The dose response curve in this International Standard parallels the Quantikine™ standard curve. To convert sample values obtained with the Quantikine™ IVD Human Epo kit to approximate NIBSC/WHO 11/170 International units, use the equation below.

Benchtop: NIBSC/WHO (11/170) approximate value (mIU/mL) = $1.3746 \times \text{Quantikine™ Human Epo value (mIU/mL)}$

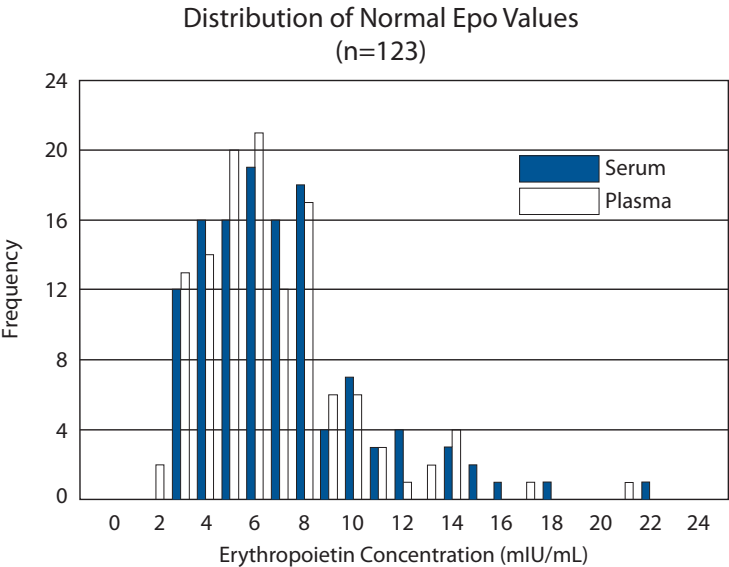
Shaker: NIBSC/WHO (11/170) approximate value (mIU/mL) = $1.4022 \times \text{Quantikine™ Human Epo value (mIU/mL)}$

Based on data generated in August 2025.

EXPECTED VALUES

Erythropoietin concentrations were obtained from 123 normal individuals from the Minneapolis/St. Paul, Minnesota area. Using the nonparametric method for the analysis of reference values outlined in the NCCLS publication *“How to Define, Determine, and Utilize Reference Intervals in the Clinical Laboratory”* (NCCLS Document C28-P; Vol. 12, No. 2) the following reference ranges (2.5-97.5 percentile) were established for Epo in serum and EDTA plasma. Each testing laboratory should establish its own normal range.

Epo Normal Ranges	
Serum	EDTA plasma
3.3-16.6 mIU/mL	3.1-14.9 mIU/mL



Patients suffering from polycythemia rubra vera may have Epo concentrations within the normal range, whereas those suffering from secondary polycythemia may have elevated concentrations of serum Epo (17). Polycythemia rubra vera patients who undergo phlebotomy may have elevated serum Epo concentrations.

Patients suffering from most anemias will present with higher than normal concentrations of serum Epo, whereas those suffering from anemia associated with chronic renal failure may have serum Epo concentrations within the normal range of this assay (18). Anemic patients who receive transfusions may exhibit lower than expected serum Epo concentrations.

Abnormally high concentrations of serum Epo may also be observed in various other pathological states including renal neoplasms, benign tumors, polycystic kidney disease, renal cysts and hydronephrosis (19).

The results of this assay should be used in conjunction with information available from clinical evaluations and other diagnostic procedures.

PERFORMANCE CHARACTERISTICS

Sensitivity

The sensitivity of the Quantikine™ IVD Human Epo ELISA minimum detectable dose is typically less than 0.6 mIU/mL. This was determined by adding two standard deviations to the mean O.D. of twenty replicates of the zero standard and calculating the corresponding concentration from the standard curve.

PERFORMANCE CHARACTERISTICS *CONTINUED*

Specificity

This assay recognizes natural and recombinant human Epo.

The factors listed below were prepared at 50 ng/mL in Specimen Diluent and assayed for cross-reactivity. Preparations of the following factors at 50 ng/mL in a mid-range Epo control were assayed for interference. No cross-reactivity or interference was observed using both the benchtop and shaker protocols.

Recombinant human:

MAN-1A1	DSC-3
CD-26	Clusterin
3-OST1	LDLR
CSPG-4	SALM-4
BMP-4	Eph B4

Recombinant human Epo Receptor did not cross-react but did interfere at concentrations > 400 pg/mL in the Quantikine™ IVD Human Epo ELISA using both the shaker and benchtop methods.

Interference of matrix-associated factors was tested by spiking a mid-range Epo control and Specimen Diluent with selected amounts of the following substances and analyzing for the presence of Epo in the Quantikine™ IVD Human Epo ELISA using both the benchtop and shaker protocols. No interference was observed.

Protein	Amount Added
Albumin	3 g/dL
Bilirubin	20 mg/dL
Gamma globulin	5 mg/dL
Hemoglobin	30 mg/dL
Transferrin	200 mg/dL
Triolein	500 mg/dL
α1-Acid Glycoprotein	80 mg/dL
α-2 Macroglobulin	30 mg/dL

PERFORMANCE CHARACTERISTICS *CONTINUED*

Accuracy

1. Recovery was estimated by addition of recombinant human Epo at three different levels throughout the range of the assay into six plasma and six serum specimens. The percent recovery of the added Epo was calculated from the equation:

$$\% \text{ recovery} = \frac{\text{measured value after addition} - \text{measured value before addition}}{\text{measured value of the added material}}$$

Mean recoveries are shown in the following table.

The recovery of human Epo spiked to levels throughout the range of the assay in various matrices was evaluated.

Specimen Type	Benchtop Protocol		Shaker Protocol	
	Average % Recovery	Range	Average % Recovery	Range
Serum (n=6)	100	84-115	101	96-107
EDTA plasma (n=6)	99	81-114	96	84-105

2. To determine whether the Quantikine™ IVD Human Epo ELISA exhibits a high dose hook, a specimen containing Epo at a concentration twenty-fold higher than the highest kit standard was serially diluted and measured in duplicate using both the benchtop and shaker assay protocols. No high dose hook was observed at 4000 mIU/mL for either protocol.

3. Six matched serum and plasma specimens samples spiked with high concentrations of human Epo were diluted with Specimen Diluent and assayed using the Quantikine™ IVD Human Epo ELISA. Diluted specimens demonstrated very good linearity when compared to neat concentrations of Epo.

Linearity

		Benchtop Protocol		Shaker Protocol	
		Serum	EDTA Plasma	Serum	EDTA Plasma
1:2	Average % of Expected	111	110	105	109
	Range (%)	106-118	104-118	99-109	101-117
1:4	Average % of Expected	109	106	102	105
	Range (%)	100-123	94-122	94-113	95-119
1:8	Average % of Expected	105	99	96	98
	Range (%)	93-123	92-118	89-106	89-114

Human Serum versus EDTA Plasma

Sixty matched serum and EDTA plasma specimens were assayed using the Quantikine™ IVD Human Epo ELISA. Results from the assay demonstrated a high correlation for benchtop protocol (r=0.9983) and for shaker protocol (r=0.9993) between the EDTA plasma and serum specimens.

PERFORMANCE CHARACTERISTICS *CONTINUED*

Assay Precision

Assay precision was assessed as outlined in the CLSI publication CLSI EP05-A3, Evaluation of Precision of Quantitative Measurement Procedures; Third Edition. The following modifications were made to the basic design described in the CLSI guideline. Three reagents lots were used instead of one. For within laboratory precision, 10 assays were run by 5 operators for each reagent lot over a period of at least 20 days with no more than 2 assays run per day.

For within assay precision, 20 replicates of each control were run for each reagent lot. Three control pools were run in both the benchtop and shaker protocols of the Quantikine™ IVD Human Epo ELISA for both within laboratory and within assay precision.

The results of each assay were tabulated and the within assay precision and within laboratory precision of each control pool was calculated.

	Benchtop Protocol						Shaker Protocol					
	Within Laboratory Precision			Within Assay Precision			Within Laboratory Precision			Within Assay Precision		
Control	1	2	3	1	2	3	1	2	3	1	2	3
Mean (mIU/mL)	9.42	40.1	116	10.6	41.3	115	9.79	40.9	116	8.81	39.4	114
n=	30	30	30	20	20	20	30	30	30	20	20	20
%CV	11.0	7.2	5.4	4.7	4.9	3.4	12.0	5.9	5.2	4.1	2.5	3.2

TROUBLESHOOTING GUIDE

Generally, assay failure is due to technical error, equipment failure or reagent failure. When an assay fails, check the expiration dates of the individual reagents and ensure that all the reagents have been stored as indicated in the product labeling. In addition, see Indications of Instability or Deterioration section for additional information. If assay performance is questionable or a problem occurs when running the assay, you may be able to isolate the problem by referring to the following table.

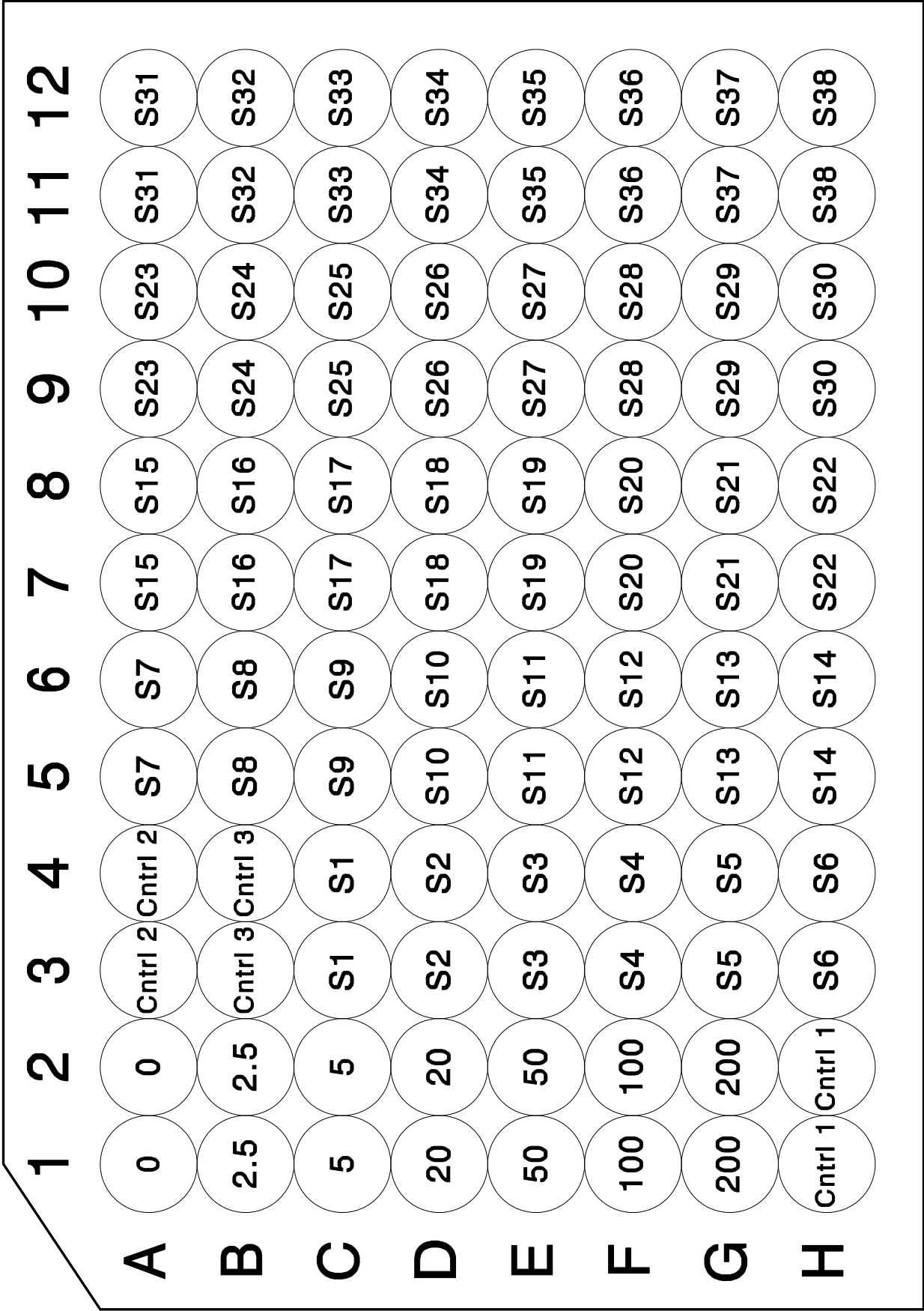
PROBLEM	POSSIBLE SOURCE	TEST OR ACTION
High % C.V.s (high variability of duplicates compared to individual laboratory precision requirements)	Incomplete washing of wells	Ensure that the wash station is working properly
	Inadequate aspiration of the wells	Wells should appear dry after aspiration
	Shaker splashing well contents onto plate cover	Calibrate the shaker to 500 ± 50 rpm
	Unequal volumes added to the wells	Ensure that the pipette is calibrated and working properly
Reduced low delta (< 0.015 O.D.) or high background	Incomplete washing of wells	Ensure that the wash station is working properly
	Inadequate aspiration of the wells	Wells should appear dry after aspiration
	Unequal volumes added to the wells	Ensure that the pipette is calibrated and working properly
Poor correlation of the Standard Curve ($r < 0.95$)	Pipetting error	Consider editing data according to individual laboratory procedures
Inadequate color development	Inadequate aspiration of the wells	Wells should appear dry after aspiration
	Unequal volumes added to the wells	Ensure that the pipette is calibrated and working properly
	Incorrect incubation times or temperatures	Adhere to recommended incubation periods and temperatures
	Conjugate or Color Reagent failure	Mix equal volumes (i.e. 100 μ L each) Epo TMB Substrate and Epo Conjugate. Blue color should develop immediately
Splashing of well contents onto adhesive plate cover	Plate shaker rpm too fast	Calibrate shaker to 500 ± 50 rpm

REFERENCES

1. Jelkmann, W. (1992) *Erythropoietin: structure, control of production, and function* in *Physiol. Reviews* **72**:449.
2. Goldwasser, E. et al. (1990) *Erythropoietin: The primary regulator of red cell formation* in **Peptide Growth Factors and their Receptors I**, Sporn, M.B. and A.B. Roberts eds., Springer-Verlag, Berlin, pp. 747.
3. Hammond, D. and S. Winnick (1974) *Paraneoplastic erythrocytosis and ectopic erythropoietins* in *Ann. N.Y. Acad. Sci.* **230**:219.
4. Chandra, M. et al. (1985) *Serum immunoreactive erythropoietin levels in patients with polycystic kidney disease as compared with other hemodialysis patients* in *Nephron* **39**:26.
5. Kenny, G.M. et al. (1970) *Erythropoietin levels in Wilms tumor patients* in *J. Urol.* **104**:758.
6. Kew, M.C. et al. (1986) *Serum erythropoietin concentrations in patients with Hepatocellular Carcinoma* in *Cancer* **58**:2485.
7. Jeffreys, R.V. et al. (1982) *Erythropoietin levels in posterior fossa haemangioblastoma* in *J. Neurol. Neurosurg. Psychiatry* **45**:264.
8. Bohling, T. et al. (1987) *Erythropoietin in capillary hemangioblastoma* in *Acta Neuropathol.* **74**:324.
9. Waldmann, T.A. and W. Rosse (1964) *Tumors producing Erythropoiesis-Stimulating factors* in **Hemoglobin, Its Precursors and Metabolites**, Sundermann, F.W. and F.W. Sundermann Jr. Eds., Lippincott, Philadelphia, pp. 276 - 280.
10. Fried, W. et al. (1968) *Leiomyoma and Erythrocytosis: A tumor producing a factor which increases Erythropoietin production. A report of case* in *Blood* **31**:813.
11. Sherwood, J.B. and E. Goldwasser (1979) *A radioimmunoassay for erythropoietin* in *Blood* **53**:885.
12. Baer, A.N. et al. (1987) *Blunted erythropoietin response to anaemia in rheumatoid arthritis* in *Br. J. Haematol.* **66**:559.
13. Spivak, J.L. et al. (1989) *Serum immunoreactive erythropoietin in HIV-infected patients* in *J. Am. Med. Assoc.* **261**:3104.
14. Miller, C.B. et al. (1990) *Decreased erythropoietin response in patients with the anemia of cancer* in *New Engl. J. Med.* **322**:1689.
15. Jelkmann, W. et al. (1990) *Modulation of the production of erythropoietin by cytokines: In vitro studies and their clinical applications* in *Contr. Nephrol.* **87**:68.
16. Cotes, P.M. and D.R. Bangham (1961) *Bio-assay of erythropoietin in mice made polycythaemic by exposure to air at a reduced pressure* in *Nature* **191**:1065.
17. Koeffler, H.P. and E. Goldwasser (1981) *Erythropoietin radioimmunoassay in evaluating patients with polycythemia* in *Ann. Intern. Med.* **94**:44.
18. Merck Manual of Diagnosis and Therapy, 14th Edition, R. Berkow, Editor.
19. Thorling, E.B. (1982) *Paraneoplastic erythrocytosis and inappropriate erythropoietin production* in *Scand. J. Haematol., Suppl.* **17**:1.

PLATE LAYOUT

A sample diagram for standards, controls, and specimens is shown below.



All trademarks and registered trademarks are the property of their respective owners.

bio•techne®

R&D Systems™ Novus Biologicals™ Tocris Bioscience™
ProteinSimple™ ACD™ Asuragen®

Global info@bio-techne.com bio-techne.com/find-us/distributors
North America TEL 800 343 7475