

	CATALOG	PRODUCT DESCRIPTION, STORAGE, AND USE
LUMINEX	LYX010	Black Microplates (96 well) (Part # 641385) - A 25 pack of black medium-binding, flat-bottom 96 well polystyrene microplates.
	LYX010/384	Black Microplates (384 well) (Part # 993671) - A 25 pack of black medium-binding, flat-bottom 384-well polystyrene microplates.
DUOSET / ELISA / ELLA	WA126	<p>Wash Buffer Concentrate (Part # 895126) - 5 vials (100 mL/vial) of a 25X concentrated solution of buffered surfactant with preservative. <i>May turn yellow over time.</i></p> <ul style="list-style-type: none"> • Store unopened at 2-8 °C. Stable for up to 30 days after opening when stored at 2-8 °C. If crystals have formed in the concentrate, warm to room temperature and mix gently until the crystals have gone into solution. • Dilute 20 mL of Wash Buffer Concentrate to a final volume of 500 mL in deionized water. • Diluted Wash Buffer may be used for up to 30 days when stored at 2-8 °C. <p>Note: Wash Buffer Concentrate contains a preservative which may cause an allergic reaction. Avoid breathing mist. Wear protective gloves, clothing, eye, and face protection. Wash hands thoroughly after handling. Refer to the SDS on our website prior to use.</p>
	DY004	<p>Reagent Diluent Concentrate 3 (5X) (Part # 895956) - 10 vials (21 mL/vial) of a 25% Tween™ 20 solution in PBS.</p> <ul style="list-style-type: none"> • Store unopened at 2-8 °C. Stable for up to 30 days after opening when stored at 2-8 °C. Make all solutions fresh or filter sterilize and store for up to 1 week at 2-8 °C. • Dilute 1:5 in PBS before use. • The reagent diluent will require additional components for some DuoSets. Refer to the product-specific package insert for more information.
	DY005	<p>Reagent Additive 1 (Part # 843566) - 1 vial (21 mL) of diluted, heat-inactivated normal goat serum.</p> <ul style="list-style-type: none"> • Store unopened at 2-8 °C. Stable for up to 30 days after opening when stored at 2-8 °C. • Add 250 µL of Reagent Additive 1 per 10 mL of 1X detection antibody or reagent diluent.
	DY006	<p>ELISA Plate-Coating Buffer (Part # 896009) - 1 vial (250 mL) of sterile-filtered 1X PBS for use as a plate-coating buffer in ELISA development.</p> <ul style="list-style-type: none"> • Contains no preservatives. • Store unopened at room temperature. Store opened at 2-8 °C. Do not use past label expiration.
	DY010	<p>Sample Activation Kit 1 (Part # 993078) - 3 vials of 1N HCl (Part # 895094, 10 mL/vial) and 3 vials of 1.2 N NaOH/0.5M HEPES (Part # 895096, 10 mL/vial). Contains no preservatives.</p> <ul style="list-style-type: none"> • Store unopened at room temperature. Do not use past kit expiration date. Stable for up to 30 days at room temperature after opening provided this is within the expiration date of the kit. • Follow activation procedure provided with each application. <p>Note: May cause skin, eye, and respiratory irritation. Avoid breathing fumes. Wear protective gloves, clothing, eye, and face protection. Wash hands thoroughly after handling. Refer to the SDS on our website prior to use.</p>
	DY990	Clear Strip-well Microplates (Part # 991427) - A 25 pack of clear high-binding, flat-bottom polystyrene microplates (360 µL/well). Each microplate consists of twelve removable strips of 8 wells and a frame.
	DY990/384	Clear Microplates (Part #608616) - A 25 pack of flat-bottom 384-well polystyrene microplates.

CATALOG	PRODUCT DESCRIPTION, STORAGE AND USE
DY991	Black Strip-well Microplates (96 well) (Part # 991426) - A 25 pack of black high-binding, flat-bottom polystyrene microplates (360 µL/well) for use in chemiluminescent assays. Each microplate consists of twelve removable strips of 8 wells and a frame.
DY992	ELISA Plate Sealers (Part # 991423) - 100 adhesive strips.
DY993	Glo Substrate Reagent Pack - 5 vials of Glo Reagent A (stabilized enhanced luminol, Part # 895868, 4 mL/vial) and 5 vials of Glo Reagent B (stabilized hydrogen peroxide, Part # 895869, 8 mL/vial). <ul style="list-style-type: none"> • Store unopened at 2-8 °C. Stable for up to 30 days after opening when stored at 2-8 °C. Use within 15 minutes to 4 hours of mixing. • This reagent may be substituted for H₂O₂/tetramethylbenzidine as a substrate for horseradish peroxidase in ELISAs. Use black microplates for chemiluminescent assays (Catalog # DY991). Mix 1 part Glo Reagent A to 2 parts Glo Reagent B in a plastic container, cap, and protect from light. • Add 100 µL of the mixture to each well and incubate at room temperature while protecting from light. Read the microplate on a luminometer within 5-20 minutes.
DY994	Stop Solution (Part # 895032) - 15 vials (6 mL/vial) of 2N sulfuric acid. <ul style="list-style-type: none"> • Store unopened at 2-8 °C. Stable for up to 30 days after opening when stored at 2-8 °C. • Add 50 µL/well to stop the action of horseradish peroxidase on the substrate, tetramethylbenzidine. If the color in the wells is green or the color change does not appear to be uniform, gently tap the plate to ensure thorough mixing. <p>Note: Wear protective gloves, clothing, eye, and face protection. Wash hands thoroughly after handling. Refer to the SDS on our website prior to use.</p>
DY994B	ELISA Stop Solution (Part # 642827) - 16 vials (12 mL/vial) of methanesulfonic acid <ul style="list-style-type: none"> • Store unopened at 2-8 °C. Stable for up to 30 days after opening when stored at 2-8 °C. • Add 100 µL/well to stop the action of horseradish peroxidase on the substrate, tetramethylbenzidine. If the color in the wells is green or the color change does not appear to be uniform, gently tap the plate to ensure thorough mixing. <p>Note: Wear protective gloves, clothing, eye, and face protection. Wash hands thoroughly after handling. Refer to the SDS on our website prior to use.</p>
DY994B-250	ELISA Stop Solution, 250 mL (Part # 642953) - 1 vial (250 mL/vial) of methanesulfonic acid <ul style="list-style-type: none"> • Store unopened at 2-8 °C. Stable for up to 30 days after opening when stored at 2-8 °C. • Add 100 µL/well to stop the action of horseradish peroxidase on the substrate, tetramethylbenzidine. If the color in the wells is green or the color change does not appear to be uniform, gently tap the plate to ensure thorough mixing. <p>Note: Wear protective gloves, clothing, eye, and face protection. Wash hands thoroughly after handling. Refer to the SDS on our website prior to use.</p>
DY995	Reagent Diluent Concentrate 2 (10X) (Part # 841380) - 5 vials (21 mL/vial) of a 10% BSA solution. <ul style="list-style-type: none"> • Store unopened at 2-8 °C. Stable for up to 30 days after opening when stored at 2-8 °C. Make all solutions fresh or filter sterilize and store for up to 1 week at 2-8 °C. For assays that require 1% BSA in PBS, dilute 1:10 in deionized water before use. • For assays that require 0.1% BSA in TBS, dilute 1:100 with Tris buffer + 0.05% Tween 20, pH 7.2-7.4. • The reagent diluent will require additional components for some DuoSets™. Refer to the product-specific package insert for more information.
DY997	Reagent Diluent Concentrate 1 (Part # 840149) - 1 vial (21 mL) of a buffered protein base. <ul style="list-style-type: none"> • Store unopened at 2-8 °C. Stable for up to 60 days after opening when stored at 2-8 °C. • Dilute 1.4 mL to a final volume of 100 mL with PBS + 0.05% Tween 20, pH 7.2-7.4. • Use the diluted reagent within 8 hours.

CATALOG	PRODUCT DESCRIPTION, STORAGE AND USE
DY998	<p>Streptavidin-HRP (Part # 890803) - 5 vials (1 mL/vial) of streptavidin conjugated to horseradish peroxidase.</p> <ul style="list-style-type: none"> • Store unopened at 2-8 °C. Stable for up to 6 months after opening when stored at 2-8 °C. Do Not Freeze. • For use in ELISAs, dilute to the working concentration specified on the vial label using 1% BSA/PBS or Reagent Diluent Concentrate 2 (Catalog # DY995) diluted 1:10 in deionized water. • For use in Western Blots, the suggested dilution is 1:200-1:2000 in an appropriate diluent. • Use the diluted reagent within 60 minutes.
DY999	<p>Substrate Reagent Pack - 8 vials of Color Reagent A (stabilized hydrogen peroxide, Part # 895000) and 8 vials of Color Reagent B (stabilized tetramethylbenzidine, Part # 895001); 12 mL/vial.</p> <ul style="list-style-type: none"> • Store unopened at 2-8 °C. Stable for up to 30 days after opening when stored at 2-8 °C. • Warm to room temperature before use. Color Reagent A and Color Reagent B should be mixed together in equal volumes within 15 minutes of use. Protect from light. Add 100 µL of the mixture to each well and incubate at room temperature for 20 minutes while protecting from light. Add Stop Solution (Catalog # DY994) and read the microplate. <p>Note: <i>Color Reagent B may cause skin, eye, and respiratory irritation. Avoid breathing fumes. Wear protective gloves, clothing, eye, and face protection. Wash hands thoroughly after handling. Refer to the SDS on our website prior to use.</i></p>
DY999B	<p>Substrate Reagent Pack - 16 vials of TMB ELISA Substrate (Part # 642736); 12 mL/vial.</p> <ul style="list-style-type: none"> • Store unopened at 2-8 °C. Stable for up to 30 days after opening when stored at 2-8 °C. • Warm to room temperature before use. Protect from light. Add 100 µL to each well and incubate at room temperature for 20 minutes while protecting from light. Add Stop Solution (Catalog # DY994B or DY994B-250) and read the microplate.
DY999B-250	<p>TMB ELISA Substrate, 250 mL - 1 bottle of TMB ELISA Substrate (Part # 642954); 250 mL/bottle.</p> <ul style="list-style-type: none"> • Store unopened at 2-8 °C. Stable for up to 30 days after opening when stored at 2-8 °C. • Warm to room temperature before use. Protect from light. Add 100 µL to each well and incubate at room temperature for 20 minutes while protecting from light. Add Stop Solution (Catalog # DY994B or DY994B-250) and read the microplate.
DYTMBS0	<p>TMB ELISA Substrate and ELISA Stop Solution Pack - 1 bottle of TMB ELISA Substrate (Part # 642954); 250 mL/ bottle and 1 bottle of ELISA Stop Solution (Part # 642953); 250 mL/ bottle.</p> <ul style="list-style-type: none"> • Store unopened at 2-8 °C. Stable for up to 30 days after opening when stored at 2-8 °C. • Warm to room temperature before use. Protect from light. Add 100 µL of TMB ELISA Substrate to each well and incubate at room temperature for 20 minutes while protecting from light. Add 100 µL/well of ELISA Stop Solution. If the color in the wells is green or the color change does not appear to be uniform, gently tap the plate to ensure thorough mixing. Measure the optical density of each well within 30 minutes using a microplate reader at 450 nm. <p>Note: <i>Wear protective gloves, clothing, eye, and face protection. Wash hands thoroughly after handling. Refer to the SDS on our website prior to use.</i></p>

DYC001

Sample Diluent Concentrate 1 (5X) (Part # 895562) - 5 vials (21 mL/vial) of a 5-fold concentrated solution containing 5X PBS, 5 mM EDTA, and 2.5% Triton X-100.

- Store unopened at 2-8 °C. Stable for up to 30 days after opening when stored at 2-8 °C. Make all solutions fresh or filter sterilize and store for up to 1 week at 2-8 °C. Do not use beyond the expiration date.
- Sample Diluent Concentrate 1 (5X) is the base buffer for the following IC Diluents and Lysis Buffers. Refer to Table 1 for detailed preparation information.

TABLE 1

Diluent Name	Sample Diluent Concentrate 1 (5X) Required	Urea Required	0.5 M NaF Required	1 M NaCl Required	Protease & Phosphatase Inhibitors Recommended*	Adjust with DI H ₂ O to Final Volume
IC Diluent #3	10 mL	3.003 g	0.5 mL	NA	No	50 mL
IC Diluent #4	10 mL	NA	NA	NA	No	50 mL
IC Diluent #7	10 mL	18.02 g	0.5 mL	NA	No	50 mL
IC Diluent #8	10 mL	NA	0.5 mL	NA	No	50 mL
IC Diluent #9	10 mL	NA	NA	25 mL	No	50 mL
Lysis Buffer #4	2 mL	NA	NA	5 mL	Yes	10 mL
Lysis Buffer #6	2 mL	3.604 g	0.1 mL	NA	Yes	10 mL
Lysis Buffer #12	2 mL	NA	NA	NA	Yes	10 mL
Lysis Buffer #13	2 mL	NA	0.2 mL	1.5 mL	Yes	10 mL

*Refer to the specific DuoSet™ IC kit insert for details.

DYC002

Sample Diluent Concentrate 2 (2X) (Part # 895891) - 1 bottle (250 mL) of a 2-fold concentrated solution containing 2% NP-40 Alternative, 40 mM Tris, 274 mM NaCl, 20% Glycerol, 4 mM EDTA, and 2 mM activated Sodium Orthovanadate. *May turn yellow over time.*

- Store unopened at 2-8 °C. Stable for up to 30 days after opening when stored at 2-8 °C. Make all solutions fresh or filter sterilize and store for up to 1 week at 2-8 °C. Do not use beyond the expiration date.
- Sample Diluent Concentrate 2 (2X) may be used as the base buffer for the following IC Diluent and Lysis Buffer. Refer to Table 2 for detailed preparation information.

TABLE 2

Diluent Name	Sample Diluent Concentrate 2 (2X) Required	25 mg/mL Leupeptin Required	Aprotinin Required	Protease Inhibitors Recommended*	Adjust with DI H ₂ O to Final Volume
IC Diluent #12	25 mL	NA	NA	No	50 mL
Lysis Buffer #9	5 mL	0.004 mL	0.1 mg	Yes	10 mL

*Refer to the specific DuoSet IC kit insert for details.