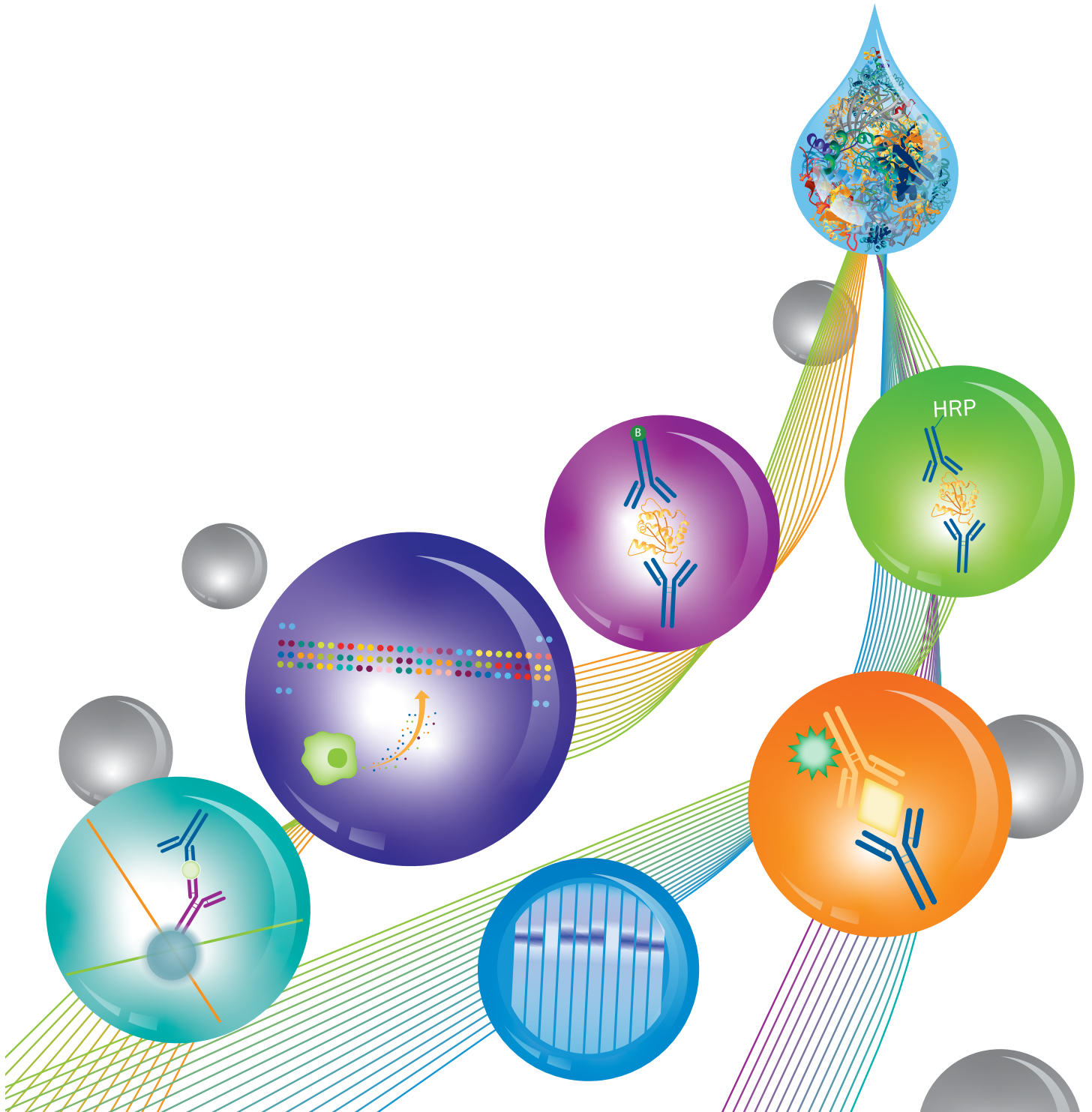


biotechne®

Quality Results Start at the Core

Bio-Techne Immunoassays — Powered by R&D Systems



What's at the Core of your Immunoassay?

Your results matter. So what's inside your immunoassay should too.

R&D Systems' antibodies and proteins are the core of every Bio-Techne immunoassay platform. Assays that are backed by over 68,000 citations and the expertise behind our gold-standard Quantikine® ELISAs. Antibodies are highly specific, manufactured in-house to ensure reproducibility and tested for suitability on every application we develop.

We've got an immunoassay platform that fits every stage of your research journey. And quality at the core means you get to accurate answers no matter what the question is. When your experimental needs change, your results won't. For most assays, R&D Systems' antibodies move with you between platforms so you can always trust your results along the way. This is done by calibrating to Quantikine kits across these immunoassays.

Screening Simplicity: Proteome Profiler Arrays

When you need: easy screening, many answers, low technology barrier



Proteome Profiler™ Antibody Arrays allow for the measurement of up to 119 proteins in a single sample. No specialized equipment required, and they are easy to use! There is a great chance you will find what you want and a whole lot more! All arrays utilize chemiluminescence for detection and select arrays are also suitable for use with the LI-COR® and FluorChem™ detection systems. Screen your samples and array the possibilities.

The DIY Solution: DuoSet ELISAs

When you need: quality results on a budget, the largest selection on the planet



Like developing your own ELISAs or need a solution that goes easy on the budget? Try our DuoSet® ELISAs. With a little DIY up front, you'll get results comparable to our other immunoassays — just prep the buffers and coat the plates. Choose from a huge menu of over 1000 single-analyte DuoSet ELISAs to measure natural, recombinant, intracellular, extracellular, or phosphorylated proteins.

The Citation Pack Leader: Quantikine ELISAs

When you need: the gold standard in single-analyte detection



Quantikine® ELISAs are the most trusted on the market, not to mention the most referenced single-analyte immunoassays to date. Quantikine assays are also master-calibrated for consistency. What does that mean for you? The results you got yesterday, last month or even last year will be the same as today's, and their 96-well assay format works with any plate reader you have in the lab. Take your pick of over 450 ready to use kits to measure a wide range of molecules including cytokines, chemokines, growth factors, proteases and more.

The Plex Builder: Luminex Assays

When you need: the most flexibility and plexibility



Plex 50 out of 450. Max out on answers for your most precious samples with our bead-based Luminex® Assays. Build a custom plex of up to 50 targets from over 450 available analytes — our Luminex Assay Customization Tool makes it simple. Here's the added bonus: you'll have access to 77 unique human analytes, 11 unique mouse analytes and 1 unique rat analyte you can't get anywhere else.

Plex by panel or analyte. Need a performance boost while you plex? Luminex High Performance Assays bump your sensitivity down to 1 pg/mL with unparalleled accuracy. Head to head testing shows that our cytokine discovery panel has superior sample recovery compared to other Luminex-based kits. Go with one of 16 preconfigured panels or build your own and multiplex up to 45 analytes. They'll let you run assays using <75 µL of sample even on complex matrices like serum, plasma, urine, saliva and milk. When you need to make a switch, Luminex High Performance Assays are correlated with our gold-standard Quantikine ELISAs so you'll get similar results across platforms.

Gel-free, blot-free, hands-free – Simple Western

When you need: fast answers, small sample size



Simple Western™ lets you separate and analyze proteins by size from 2–440 kDa either by immunoassay or total protein analysis. And there's even a charge-based assay if you're doing post-translational modification characterization. You'll get quantitative results, reproducibility that's spot on, and use less sample in the process. So what's the big difference between Simple Western and everything else? It's the only gel-free, blot-free, hands-free capillary-based immunoassay platform that integrates and automates the entire protein separation and detection process. So you'll have more time to get down to science.


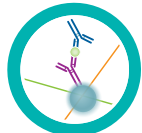




Your Next Generation ELISA: Simple Plex Assays

When you need: mega speed, max sensitivity, awesome reproducibility



Why not leave the complexity of running immunoassays to us? Automated Simple Plex™ assays on E11a let you run ELISAs hands-free. Just load the cartridge and go, answers are ready in an hour. You'll also get sub pg/mL detection and reproducibility in the low single digits — even across multiple users and labs. Run a single analyte on up to 72 samples or 4-plex it on up to 32 samples with zero cross-reactivity. And you'll only use 25 µL of sample at most. Because Simple Plex assays are correlated against our gold-standard Quantikine ELISAs, it's always easy to transfer to another platform whenever you need to.

Which Immunoassay is Right for You?

	Description	Format	Quantitative	Sample volume	Number of analytes
	Proteome Profiler™ Arrays	Membrane	Relative	200–500 µL	15–119
	Luminex® Assays	Bead	✓	25–50 µL	Up to 50
	Simple Western™	Capillary Cartridge	✓	3 µL	>1
	DuoSet® ELISAs	Flexible	✓	100 µL	1
	Quantikine® ELISAs	96-well plate	✓	10–200 µL	1
	Simple Plex™ Assays	Cartridge	✓	2.5–25 µL	Up to 4

R&D SYSTEMS

NOVUS BIOLOGICALS

TOCRIS

protein**simple**

A&C D

bio-techne

Global info@bio-techne.com bio-techne.com/find-us/distributors TEL +1 612 379 2956
 North America TEL 800 343 7475 Europe | Middle East | Africa TEL +44 (0)1235 529449
 China info.cn@bio-techne.com TEL +86 (21) 52380373

bio-techne.com

