



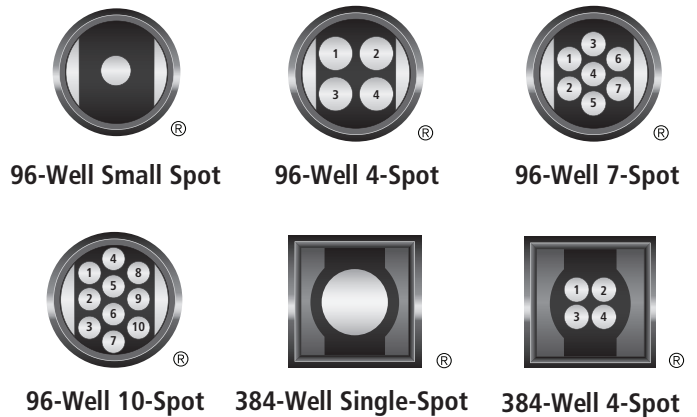
MSD[®] Prototype Printing Services

MESO SCALE DISCOVERY[®] offers Prototype Printing Services to coat MSD plates with materials of your choice (antibodies, proteins, peptides, carbohydrates, polysaccharides, or oligonucleotides). These services provide a rapid and convenient way to convert ELISAs or create multiplex assays from your own or commercially available materials. MSD GOLD[™] Prototype Printing Service includes additional characterization of antibodies, which accelerates the development of assays and helps manage variability originating from changes in raw material that can occur from lot to lot. MSD's Prototype Printing Services leverage our ISO 9001:2008-certified quality management system and extensive experience in manufacturing, providing researchers with a cost-effective and convenient route to high-quality coated plates.

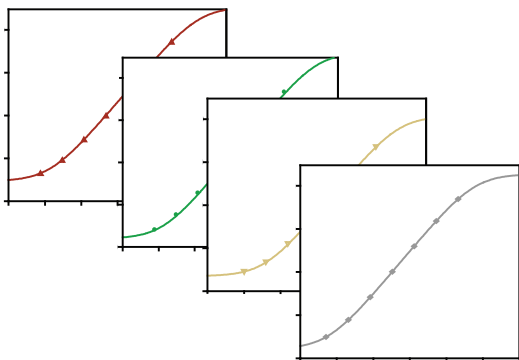
Proven Benefits of MSD Assays

- “[Compared] with the ELISA, the Pn [pneumococcal] ECL assay provides a **25-fold reduction in required serum volumes** for the reference standard and test samples and a **200-fold reduction** in the amount of Ps [polysaccharides] required.” (Marchese et al. 2009)
- “Quantitative assessment of IGF-IR in patient samples and cell lines was performed using a novel sandwich assay (ECL) as it provides **sensitivity (~1 pg), accuracy, reproducibility, and absolute quantification.**” (Cao et al. 2008)
- “The assays are **robust**...and offer **significant advantages (in terms of speed and quantitation)** over western blots.” (Gowan et al. 2007)

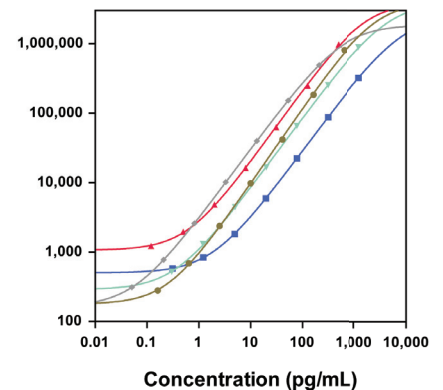
MSD Spot Patterns for Printing Services



Multiplex with MSD Prototype Printing Services



Conserve Sample
Improve Workflow
Large Dynamic Range
Exquisite Sensitivity

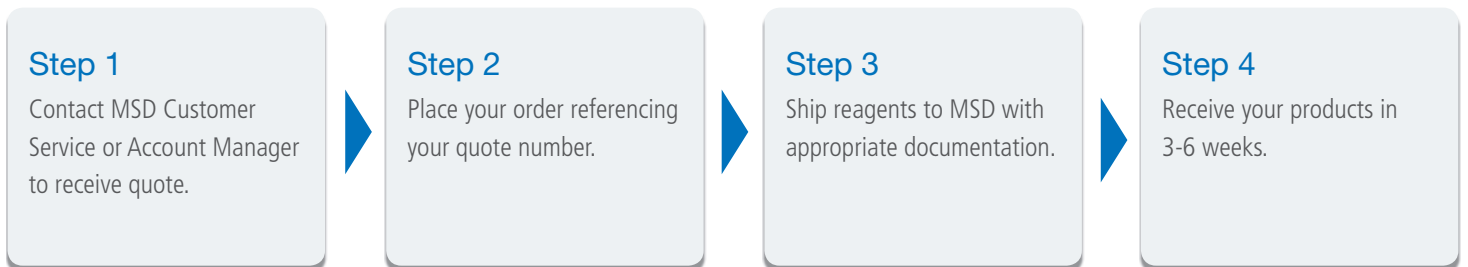


Service	MSD GOLD Prototype Printing Service	MSD Prototype Printing Service
Antibody buffer exchanged to remove interferences	✓	
Add-on size exclusion chromatography to remove aggregates (optional)	✓	
Pre-printing analytical reagent characterization	✓	
Plate printing under controlled environment	✓	✓
In-process QC including images of fluids dispensed	✓	✓
Plate packaging in single plate pouches	✓	✓
Quality control for coating uniformity	✓	✓
Certificate of Analysis	Includes characterization data	Upon request

Quality Control (QC) for Coating Uniformity

Plates are manufactured utilizing documented procedures controlled under MSD's Quality Management System (QMS) using precise, low volume dispensers and automation systems. MSD's QMS has received ISO 9001:2008 certification for the design, development, manufacture, distribution and service of instruments, plates, reagents, assays and kits for research purposes. MSD has stringent QC processes built into various stages of plate printing, including testing for electrical conductivity in every well, visual inspection during dispensing, and testing for antibody coating consistency.

Contact us to discuss your needs.



Phone: 1-240-314-2795 . Fax: 1-301-990-2776
Email: customerservice@mesoscale.com

MESO SCALE DISCOVERY, MSD, www.mesoscale.com, MSD GOLD, SMALL SPOT (design), 96 WELL 1, 4, 7, & 10-SPOT (designs), 384 WELL 1 & 4-SPOT (designs), MSD (design), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.
 © 2015 Meso Scale Diagnostics, LLC. All rights reserved.



For Research Use Only. Not for use in diagnostic procedures.
 MK-DS-220-v1-2015Jan