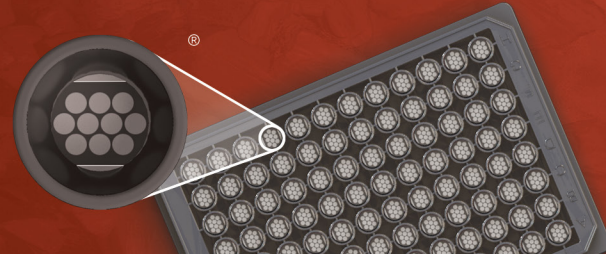


# MSD<sup>®</sup> Proinflammatory Panel 1 (rat) Control Pack



## Ordering Information

MSD Customer Service  
Phone: 1-240-314-2795  
Fax: 1-301-990-2776  
Email: CustomerService@mesoscale.com  
[www.mesoscale.com/support](http://www.mesoscale.com/support)

## Scientific Support

Phone: 1-240-314-2798  
Email: ScientificSupport@mesoscale.com  
[www.mesoscale.com/support](http://www.mesoscale.com/support)

## Company Address

MESO SCALE DISCOVERY<sup>®</sup>  
A division of  
Meso Scale Diagnostics, LLC.  
1601 Research Boulevard  
Rockville, MD 20850-3173 USA

[www.mesoscale.com](http://www.mesoscale.com)<sup>®</sup>

**For Research Use Only.  
Not for use in diagnostic  
procedures.**

<b>Catalog #</b>	C4044-1
<b>Contents</b>	Proinflammatory Panel 1 (rat) Control 1 (5 vials,* lyophilized) Proinflammatory Panel 1 (rat) Control 2 (5 vials,* lyophilized) Proinflammatory Panel 1 (rat) Control 3 (5 vials,* lyophilized) *Supplied as part of a V-PLEX <sup>®</sup> Plus kit, quantity will match number of plates ordered.
<b>Storage</b>	2–8°C

## Summary and Intended Use

The use of controls to monitor the analytical performance and reliability of test methods is an essential component of good laboratory practice. The Proinflammatory Panel 1 (rat) Control Pack consists of 3 levels of controls, each containing known concentrations of all rat cytokines that are detected by the V-PLEX Proinflammatory Panel 2 (rat) Kit. Proinflammatory Panel 1 (rat) Controls 1, 2, and 3 are prepared by spiking recombinant calibrators into a non-rat matrix diluent. The controls are supplied as lyophilized powder.

## Storage and Handling

To maximize the consistency of measured values across vials, the controls must be stored at the temperature recommended above. These controls are a one-time use product; discard any unused reconstituted product.

To use, first reconstitute each vial of Proinflammatory Panel 1 (rat) control in 250  $\mu$ L of Diluent 42. Do not mix or vortex the vials. Wait for a minimum of 15-30 minutes, briefly vortex, and then dilute to match the sample dilution (4-fold is the recommended sample dilution for this kit). Add diluted control solutions directly to the MSD Proinflammatory Panel 2 (rat) plate and assay as unknown samples. Discard unused diluted control material.

## Safety

Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling controls. Handle and dispose of all hazardous samples properly in accordance with local, state, and federal guidelines. Additional product-specific safety information is available in the safety data sheet, which can be obtained from MSD Customer Service.

## Assignment of Control Values

The controls are provided to assess reproducibility of assay performance, and precision CVs are expected to be less than 25%. The certificate of analysis contains the concentrations of the controls measured at MSD across three lots. Even with good laboratory practices, site-to-site differences may occur; therefore, to establish accuracy specifications, it is recommended that each lab establish its own nominal values and acceptance range for the control concentrations.

MESO SCALE DISCOVERY, Meso Scale Diagnostics, MSD, V-PLEX, www.mesoscale.com, MSD (design), 96 WELL 10-SPOT (design), and Spot the Difference are trademarks and/or service marks of Meso Scale Diagnostics, LLC.  
©2013-2014, 2022 Meso Scale Diagnostics, LLC. All rights reserved.