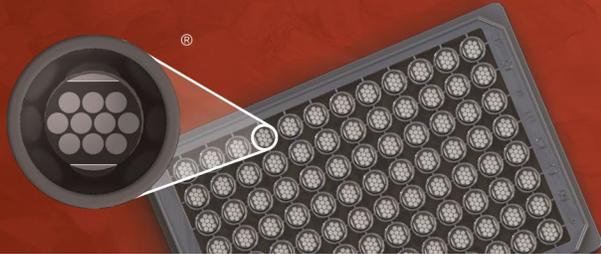


Vascular Injury Control Pack 1



www.mesoscale.com[®]

Ordering Information

MSD Customer Service
Phone: 1-240-314-2795
Fax: 1-301-990-2776
Email: CustomerService@mesoscale.com

www.mesoscale.com/support

Scientific Support

Phone: 1-240-314-2798
Email: ScientificSupport@mesoscale.com

www.mesoscale.com/support

Company Address

MESO SCALE DISCOVERY[®]
A division of
Meso Scale Diagnostics, LLC.
1601 Research Boulevard
Rockville, MD 20850-3173 USA

Catalog No	C4687-1
Contents	Vascular Injury Control 1 (5 vials,* 250 µL/vial) Vascular Injury Control 2 (5 vials,* 250 µL/vial) Vascular Injury Control 3 (5 vials,* 250 µL/vial) *Supplied as part of a V-PLEX [®] Plus kit, quantities will match the number of plates ordered.
Storage	≤ -70°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 Vascular Injury Control Pack 1 consists of 3 levels of controls, each containing known concentrations of all analytes that are detected by the V-PLEX Plus Vascular Injury Panel 2 (human) Kit (catalog #K15198G-1). The controls span the linear ranges of the assays in the V-PLEX Plus Vascular Injury Panel 2 (human) Kit and are designed to test the analytical range of the assays. Vascular Injury Controls 1, 2, and 3 are prepared by spiking recombinant proteins into diluent. The controls are supplied as frozen liquids.

Note: The control pack containing native CRP protein (catalog number: C4198-1) was discontinued and replaced by a control pack containing recombinant CRP protein (catalog number: C4687-1).

Storage and Handling

To maximize the consistency of measured values across vials, the controls must be stored at the temperature recommended above. The controls can tolerate 3 freeze-thaw cycles. Discard unused material after the third freeze-thaw cycle. Controls do not require dilution before use. Add a neat solution directly to the MSD Vascular Injury Panel 2 (human) plate and assay as unknown samples.

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 the 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 should establish its own nominal values and acceptance range for the control concentrations.

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

MESO SCALE DISCOVERY, Meso Scale Diagnostics, MSD, WWW.MESOSCALE.COM, MSD (design), V-PLEX, 96 WELL 10-SPOT (design), and Spot the Difference are trademarks and/or service marks of Meso Scale Diagnostics, LLC.
©2023 Meso Scale Diagnostics, LLC. All rights reserved.



MK-DS-1165-v1-2023Feb