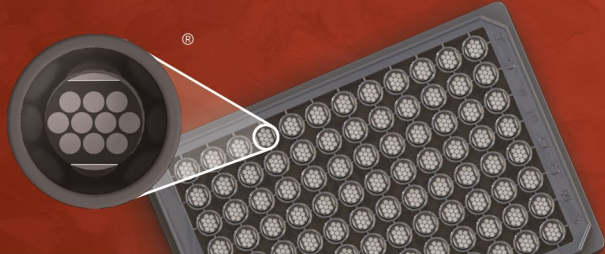


Vascular Injury Panel 2 (human) Calibrator Blend



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	C0687-2
Contents	Vascular Injury Panel 2 (human) Calibrator Blend (1 vial; 20 µL)
Storage	≤70°C

Summary and Intended Use

The Vascular Injury Panel 2 (human) Calibrator Blend is a frozen liquid combination of 4 proteins: SAA, CRP, VCAM-1, and ICAM-1. Vascular Injury Panel 2 (human) Calibrator Blend was developed for use as an assay calibrator for MSD assays in the V-PLEX® and V-PLEX Plus Vascular Injury Panel 2 (human) Kits (catalog numbers: K15198D and K15198G). The utility of this product has not been evaluated in other applications.

Note: The calibrator blend containing native CRP protein (catalog number: C0198-2) was discontinued and replaced by a calibrator blend containing recombinant CRP protein (catalog number: C0687-2).

Reagent

Vascular Injury Panel 2 (human) Calibrator Blend includes recombinant proteins. SAA protein is derived from a wheat germ expression system, CRP from a mammalian cell line, VCAM-1 from Chinese hamster ovary cells, and ICAM-1 from a mouse myeloma cell line. Vascular Injury Panel 2 (human) calibrator blend is supplied as a frozen liquid that yields a top-of-curve calibrator concentration upon a 20-fold dilution into assay diluent (Diluent 101; catalog #R51AD).

Storage and Handling

To maximize consistency in calibrator values across assay runs, the calibrator must be stored at the recommended temperature and handled per the instructions provided in this document. Thaw the calibrator blend on ice before initiating serial dilutions. The material is intended for one-time use, however, the calibrator blend can tolerate 3 freeze-thaw cycles.

Vascular Injury Panel 2 (human) calibration curves are constructed based on the concentration of calibrator in each well. Refer to the assay-specific product insert for calibrator preparation instructions. Excess diluted calibrator should be discarded after use.

Safety

Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling calibrators. 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.

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-1164-v1-2023Feb