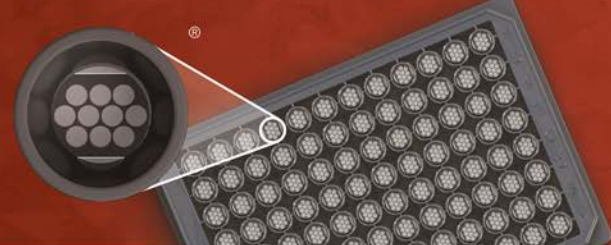


MSD® Vascular Injury Panel 2 (human)

Calibrator Blend



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

www.mesoscale.com®

Catalog No	C0198-2
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™ Vascular Injury Panel 2 (human) Kit (catalog #K15198D-1). The utility of this product has not been evaluated in other applications.

Reagent

Vascular Injury Panel 2 (human) Calibrator Blend includes 3 full-length recombinant proteins. SAA protein is derived from a wheat germ expression system, VCAM-1 from Chinese Hamster Ovary cells, and ICAM-1 from a mouse myeloma cell line. CRP native protein is isolated from human pleural/ascitic fluid or plasma. Vascular Injury Panel 2 (human) calibrator blend is supplied as a frozen liquid that yields top-of-curve calibrator concentration upon a 20-fold dilution into assay diluent (Diluent 101; catalog #R51AD).

Storage and Handling

In order 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 prior to 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

The Vascular Injury Panel 2 (human) Calibrator Blend contains native CRP protein and is a biosafety level 2 (BSL-2) product. Tissue donors have been tested and found negative for HBsAg, HIV-1 and HIV-2 antibodies, and HCV. 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.