

PRODUCT INFORMATION



Vaspin (human) Monoclonal Antibody (Clone VP63)

Item No. 10812

Overview

Contents: This vial contains a 1 mg/ml solution of antibody in PBS, pH 7.4.
Synonyms: OL-64, Serpin A12, Visceral Adipose Tissue-Derived Serine Protease Inhibitor
Immunogen: Recombinant human vaspin
Species Reactivity: (+) Human vaspin

Properties

Storage: -20°C (as supplied)
Stability: 2 years
Clone: VP63
Host: Mouse

Applications

Western blot: Suggested starting dilution is 1:2,000-1:5,000
ELISA (direct or indirect): Suggested starting dilution is 1:2,000-1:5,000

Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Images



Lane 1: Recombinant human vaspin (His-tagged)
Lane 2: Recombinant human vaspin (FLAG[®]-tagged)

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/03/2015

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM