

PRODUCT INFORMATION



Scavenger Receptor B2/CD36 Blocking Peptide

Item No. 300011

Overview and Properties

Contents:	This vial contains 200 µg of lyophilized peptide.
Synonyms:	Fatty Acid Translocase, GPIIb, GPIV, PAS IV, Platelet Collagen Receptor, Platelet Glycoprotein 4, Thrombospondin Receptor
Form:	Solid
Storage:	-20°C (as supplied)
Stability:	≥3 years

Procedures

This vial contains 200 µg of lyophilized peptide. The peptide is derived from the human CD36 sequence amino acids 98-114 (AKENVTQDAEDNTVSF). This blocking peptide can be used in conjunction with Cayman's Scavenger Receptor B2/CD36 Polyclonal Antibody (Item No. 100011) to block protein-antibody complex formation during immunochemical analysis of CD36.

To block antibody/protein complex formation, the following procedure is recommended:

1. Mix the Scavenger Receptor B2/CD36 Polyclonal Antibody (Item No. 100011) and blocking peptide together in a 1:1 (v/v) ratio in a microfuge tube. For example, mix 20 µl of antibody and 20 µl of peptide.*
2. Incubate for one hour at room temperature with occasional mixing prior to further dilution and application of the mixture to the immunoblot.
3. Dilute the mixture to the final working antibody concentration and apply to the slide or membrane as usual.

*This is a recommended mixture. The minimum amount of peptide needed for complete blocking has not been precisely determined and may vary depending on the sample being analyzed. The amount of peptide required may need to be increased if sufficient blocking does not occur.

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, 12/07/2023

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