

PRODUCT INFORMATION



Serum Retinol Binding Protein 4 Blocking Peptide

Item No. 10007682

Overview and Properties

Contents:	This vial contains 200 µg of peptide.
Synonyms:	sRBP4 Blocking Peptide, pRBP Blocking Peptide, Plasma Retinol Binding Protein 4 Blocking Peptide
Form:	Liquid
Storage:	-20°C (as supplied)
Storage Buffer:	200 µl TBS, pH 7.4, containing 0.1% BSA and 0.02% sodium azide
Stability:	≥3 years

Procedures

This vial contains 200 µg peptide in 200 µl TBS, pH 7.4, containing 0.1% BSA and 0.02% sodium azide. The sRBP blocking peptide (human amino acids 28-37) can be used in conjunction with Cayman's sRBP4 Polyclonal Antibody (Item No. 10007681) to block protein-antibody complex formation during immunochemical analysis of sRBP4.

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

1. Mix the sRBP4 Polyclonal Antibody (Item No. 10007681) 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.

Description

Serum retinol binding protein (sRBP4) binds one equivalent of vitamin A and is one of the major retinol carriers found in the blood of mammals.^{1,2} Human RBP4 is a monomeric 21 kDa β-sheet-rich protein that contains three disulfide bonds and belongs to the lipocalin protein family.³ sRBP4 is synthesized and sequestered in hepatocytes until retinol binding triggers its secretion.³ In plasma, sRBP4 typically forms a 1:1 complex with the 55 kDa tetrameric protein transthyretin (TTR) which prevents RBP from being removed from the plasma by glomerular filtration.⁴ Recent studies have shown that sRBP4 is an adipocyte-derived "signal" that may contribute to the pathogenesis of type 2 diabetes.^{5,6} Elevation of sRBP4 causes systemic insulin resistance whereas reduction of serum concentrations improves insulin action.⁵⁻⁸ The highest known concentrations of this protein exist in serum, liver, and skeletal muscle.^{5,8,9}

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, 10/17/2023

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

PRODUCT INFORMATION



References

1. Noy, N. *Biochem. J.* **348**(Pt. 3), 481-495 (2000).
2. Xie, Y., Lashuel, H.A., Miroy, G.J., et al. *Protein Expr. Purif.* **14**(1), 31-37 (1998).
3. Cowan, S.W., Newcomer, M.E., and Jones, T.A. *Proteins* **8**(1), 44-61 (1990).
4. Sivaprasadarao, A. and Findlay, J.B.C. *Biochem. J.* **296**(Pt. 1), 209-215 (1993).
5. Yang, Q., Graham, T.E., Mody, N., et al. *Nature* **436**(7049), 356-362 (2005).
6. Muoio, D.M. and Newgard, C.B. *Nature* **436**(7049), 337-338 (2005).
7. Graham, T.E., Yang, Q., Blüher, M., et al. *N. Engl. J. Med.* **354**(24), 2552-2563 (2006).
8. Polonsky, K.S. *N. Engl. J. Med.* **354**(24), 2596-2598 (2006).
9. Tamori, Y., Sakaue, H., and Kasuga, M. *Nat. Med.* **12**(1), 30-31 (2006).

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