

PRODUCT INFORMATION

Serum Retinol Binding Protein 4 (human recombinant)

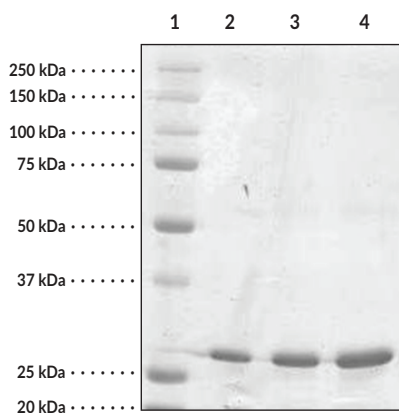
Item No. 10007818

Overview and Properties

Synonyms: RBP4, sRBP4
Source: Recombinant His-tagged protein purified from *E. coli*
Molecular Weight: ~21 kDa
Storage: -80°C (as supplied)
Stability: ≥6 months
Purity: ≥95% estimated by SDS-PAGE
Supplied in: 50 mM sodium phosphate, pH 8.2, with 100 mM sodium chloride and 20% glycerol
Protein
Concentration: *batch specific* mg/ml

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Lane 1: MW Markers
 Lane 2: 0.5 µg
 Lane 3: 1.0 µg
 Lane 4: 2.0 µg

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, 04/25/2022

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



Description

Human serum retinol binding protein 4 (sRBP4) binds to one equivalent of vitamin A and is one of the major retinol carriers found in the blood of mammals.^{1,2} Human sRBP4 is a monomeric 21 kDa β -sheet-rich protein that contains three disulfide bonds and belongs to the lipocalin protein family.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 reduced serum concentrations of sRBP4 improves insulin action.^{5,7,8} Cayman's human recombinant sRBP4 contains a C-terminal hexahistidine tag. The purified protein was characterized for its retinol binding activity. sRBP4 is stable at -20°C for six months. For long term storage aliquot the protein and store at -80°C.

References

1. Noy, N. Retinoid-binding proteins: Mediators of retinoid action. *Biochem. J.* **348**(Pt. 3), 481-495 (2000).
2. Xie, Y., Lashuel, H.A., Miroy, G.J., *et al.* Recombinant human retinol-binding protein refolding, native disulfide formation, and characterization. *Protein Expr. Purif.* **14**(1), 31-37 (1998).
3. Cowan, S.W., Newcomer, M.E., and Jones, T.A. Crystallographic refinement of human serum retinol binding protein at 2Å resolution. *Proteins* **8**(1), 44-61 (1990).
4. Sivaprasadarao, A. and Findlay, J.B.C. Expression of functional human retinol-binding protein in *Escherichia coli* using a secretion vector. *Biochem. J.* **296**(Pt. 1), 209-215 (1993).
5. Yang, Q., Graham, T.E., Mody, N., *et al.* Serum retinol binding protein 4 contributes to insulin resistance in obesity and type 2 diabetes. *Nature* **436**(7049), 356-362 (2005).
6. Muoio, D.M. and Newgard, C.B. Metabolism: A is for adipokine. *Nature* **436**(7049), 337-338 (2005).
7. Graham, T.E., Yang, Q., Blüher, M., *et al.* Retinol-binding protein 4 and insulin resistance in lean, obese, and diabetic subjects. *N. Engl. J. Med.* **354**(24), 2552-2563 (2006).
8. Polonsky, K.S. Retinol-binding protein 4, insulin resistance, and type 2 diabetes. *N. Engl. J. Med.* **354**(24), 2596-2598 (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