

PRODUCT INFORMATION



Small Integral Membrane Protein 11 (human, recombinant)

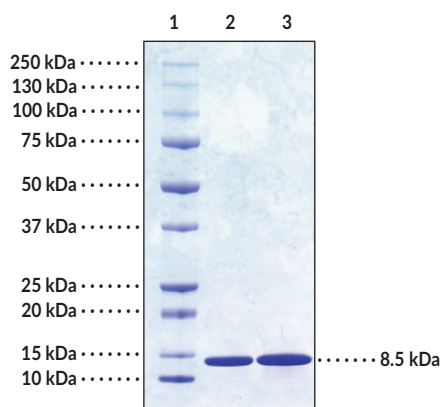
Item No. 37620

Overview and Properties

Synonyms:	Chromosome 21 Open Reading Frame 51, C21orf51, SMIM11, SMIM11A, SMIM11B, FAM165B, Family with Sequence Similarity 165, Member B
Source:	Recombinant human C-terminal Rho-1D4-tagged SMIM11 expressed in insect cells
Amino Acids:	1-58 (full length)
Uniprot No.:	P58511
Molecular Weight:	8.5 kDa
Storage:	-80°C (as supplied)
Stability:	≥6 months
Purity:	≥90% estimated by SDS-PAGE
Supplied in:	10 mM Sodium phosphate, pH 7.0, with 150 mM sodium chloride and 10% glycerol
Protein	
Concentration:	<i>batch specific</i> 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: Small Integral Membrane Protein 11 (2 µg)

Lane 3: Small Integral Membrane Protein 11 (4 µg)

SDS-PAGE Analysis of Small Integral Membrane Protein 11.

Representative gel image shown; actual purity may vary between each batch.

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, 07/06/2023

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



Description

Small integral membrane protein 11 (SMIM11) is a single-pass membrane protein encoded by the *SMIM11* gene found on human chromosome 21 (HC21), the chromosome that is trisomic in Down syndrome.¹ It is composed of a single α helical domain and is expressed in a variety of human and mouse tissues. *SMIM11A* expression is increased in an *in vitro* cancer cell model of vascular mimicry, and its silencing *via* siRNA prevents the formation of tubular structures in the same model.²

References

1. Reymond, A., Friedli, M., Henrichsen, C.N., *et al.* From PREDs and open reading frames to cDNA isolation: Revisiting the human chromosome 21 transcription map. *Genomics* **78**(1-2), 46-54 (2001).
2. Mingo, G., Pradenas, J., Babbitt, N., *et al.* SMIM11A, a novel protein involved in the mechanism of vasculogenic mimicry formation *in vitro*. *Cancer Res.* **82**(12 Suppl), 3841 (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