

PRODUCT INFORMATION



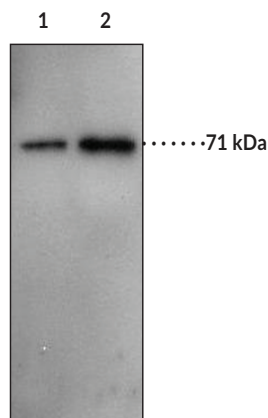
Prostaglandin Transporter (C-Term) Polyclonal Antibody

Item No. 11860

Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity purified polyclonal antibody.
Synonyms:	OATP2A1, PGT, SLCO2A1, Solute carrier organic anion transporter family member 2A1
Immunogen:	Synthetic peptide from the C-terminal region of human PGT
Species Reactivity:	(+) Human, cow, and sheep; other species not tested
Uniprot No.:	Q92959
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Applications:	Western blot (WB); the recommended starting dilution is 1:200. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Image



Lane 1: Human platelet lysate (25 µg)
Lane 2: Human platelet lysate (60 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, 11/07/2023

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

PRODUCT INFORMATION



Description

Prostaglandins (PGs) are signaling molecules that modulate physiological or pathologic functions such as platelet aggregation, vascular tone, gastric cytoprotection, uterine contraction, inflammation, cancer and Alzheimer's disease.¹⁻⁶ PG transporter (PGT) is generally expressed in endothelial and epithelial cells. Transport of extracellular PGs into cells occurs via a specific PGT, allowing PG oxidation by 15-hydroxy PG dehydrogenase or binding with their respective receptor(s).⁷⁻⁹ Cayman's peptide-affinity purified PGT polyclonal antibody detects an 85 kDa band on WB. Loss of PGT detection may occur due to cell membrane aggregation during exposure of some samples to high temperature processing.¹⁰ Treatment of the protein sample at 37°C for 15 minutes (water bath or heating block), and if desired, in parallel with another aliquot at 95°C is recommended prior to electrophoresis and blotting. When using the antibody for ICC or IF, cell permeabilization is necessary due to the intracellular location of the antibody binding site (C-terminal amino acids).

References

1. Samuelsson, B., Granström, E., Green, K., *et al.* Prostaglandins. *Annu. Rev. Biochem.* **44**, 669-695 (1975).
2. Holm, L. and Jägare, A. Role of prostaglandins in regulation of gastric mucosal blood flow and acid secretion. *Am. J. Physiol.* **263(4 Pt 1)**, G446-G451 (1992).
3. Gobbetti, A., Zerani, M., Vannucci, F., *et al.* Prostaglandins and sex steroids from corpora leutea of pregnant mares. *In vitro* studies. *Prostaglandins* **48(4)**, 262-272 (1994).
4. Harris, S.G., Padilla, J., Koumas, L., *et al.* Prostaglandins as modulators of immunity. *Trends Immunol.* **23(3)**, 144-150 (2002).
5. Engblom, D., Ek, M., Saha, S., *et al.* Prostaglandins as inflammatory messengers across the blood-brain barrier. *J. Mol. Med. (Berl)* **80(1)**, 5-15 (2002).
6. Tilley, S.L., Coffman, T.M., and Koller, B.H. Mixed messages: Modulation of inflammation and immune responses by prostaglandins and thromboxanes. *J. Clin. Invest.* **108(1)**, 15-23 (2001).
7. Kanai, N., Lu, R., Satriano, J.A., *et al.* Identification and characterization of a prostaglandin transporter. *Science* **268(5212)**, 866-869 (1995).
8. Lu, R., Kanai, N., Bao, Y., *et al.* Cloning, *in vitro* expression, and tissue distribution of a human prostaglandin transporter cDNA (hPGT). *J. Clin. Invest.* **98(5)**, 1142-1149 (1996).
9. Subbaramaiah, K., Hudis, C.A., and Dannenberg, A.J. The prostaglandin transporter regulates adipogenesis and *Aromatase* transcription. *Cancer Prev. Res. (Phila)* **4(2)**, 194-206 (2011).
10. Schägger, H. Tricine-SDS-PAGE. *Nat. Protoc.* **1(1)**, 16-22 (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