

PRODUCT INFORMATION



Prostaglandin D Synthase (hematopoietic-type; human, recombinant)

Item No. 10006593

Overview and Properties

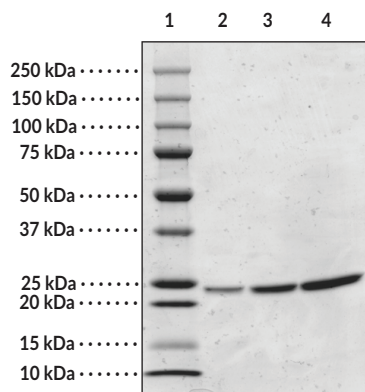
Synonyms: H-PGDS, hH-PGDS, Hematopoietic-PGDS, PGD Synthase (hematopoietic-type; human recombinant)
Source: Active recombinant human N-terminal His-tagged H-PGDS expressed in *E. coli*
Amino Acids: 1-199 (full length)
Uniprot No.: O60760
Molecular Weight: 24.3 kDa
Storage: -80°C (as supplied)
Stability: ≥2 years
Purity: *batch specific* (≥95% estimated by SDS-PAGE)
Supplied in: 50 mM sodium phosphate, pH 7.2, containing 20% glycerol, 100 mM sodium chloride, 1 mM DTT, and 0.5 mM EDTA

Protein

Concentration: *batch specific* mg/ml
Specific Activity: *batch specific* U/mg
Unit Definition: One unit of enzyme produces 1 μmole of PGD₂ per minute at 25°C in 100 mM Tris-HCl, pH 8.0, 1 mM GSH, 1 mM magnesium chloride and 40 μM PGH₂.

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

Image



Lane 1: MW Markers
Lane 2: hH-PGDS (1 μg)
Lane 3: hH-PGDS (2 μg)
Lane 4: hH-PGDS (4 μg)

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, 10/11/2021

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

Prostaglandin D synthase (PGDS) is a glutathione-dependent enzyme and member of the sigma class of glutathione-S-transferases (GSTs) that catalyzes the conversion of PGH₂ (Item No. 17020) to PGD₂ (Item No. 12010), an eicosanoid that has numerous biological functions, including vasorelaxation, recruitment of inflammatory cells, and inhibition of platelet aggregation.¹⁻³ There are two types of PGDS: lipocalin PGDS (L-PGDS; Item Nos. 10006788 | 10006787 | 10010548) and hematopoietic PGDS (H-PGDS; Item Nos. 10006593 | 10004347).³ H-PGDS is found in peripheral tissues and immune cells, including Th2 cells, antigen-presenting cells, mast cells, megakaryocytes, and eosinophils, where it is localized to the cytosol.² H-PGDS activity is increased by a variety of stimuli, including LPS, anti-IgE antibodies, phorbol 12-myristate 13-acetate (TPA; Item No. 10008014), ionomycin (Item No. 10004974), and inflammatory cytokines such as IL-13, IL-3, or IL-4.³ siRNA silencing of *Hpgds* decreases LPS-induced production of PGD₂ in mouse bone marrow-derived macrophages (BMDMs).⁴ Transgenic overexpression of *HPGDS* in mice increases croton oil-induced ear swelling and PGD₂ production, and genome-wide deletion of *Hpgds* exacerbates hypotension and vascular permeability in a mouse model of anaphylaxis.^{5,6} H-PGDS protein levels are increased in the nasal mucosa of patients with allergic rhinitis, and *HPGDS* SNPs have been found in individuals with asthma.^{1,3} Cayman's Prostaglandin D Synthase (hematopoietic-type; human, recombinant) can be used for enzyme activity assays.

References

1. Kanaoka, Y. and Urade, Y. Hematopoietic prostaglandin D synthase. *Prostaglandins Leukot. Essent. Fatty Acids* **69(2-3)**, 163-167 (2003).
2. Thurairatnam, S. Hematopoietic prostaglandin D synthase inhibitors. *Prog. Med. Chem.* **51**, 97-133 (2012).
3. Rittchen, S. and Heinemann, A. Therapeutic potential of hematopoietic prostaglandin D₂ synthase in allergic inflammation. *Cells* **8(6)**, 619 (2019).
4. Zhao, G., Yu, R., Deng, J., *et al.* Pivotal role of reactive oxygen species in differential regulation of lipopolysaccharide-induced prostaglandins production in macrophages. *Mol. Pharmacol.* **83(1)**, 167-178 (2013).
5. Sarashina, H., Tsubosaka, Y., Omori, K., *et al.* Opposing immunomodulatory roles of prostaglandin D₂ during the progression of skin inflammation. *J. Immunol.* **192(1)**, 459-465 (2014).
6. Nakamura, T., Fujiwara, Y., Yamada, R., *et al.* Mast cell-derived prostaglandin D₂ attenuates anaphylactic reactions in mice. *J. Allergy Clin. Immunol.* **140(2)**, 630-632 (2017).

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