## **PRODUCT** INFORMATION



Prostaglandin D Synthase (lipocalin-type) Polyclonal Antibody Item No. 160003

#### **Overview and Properties**

Contents:	This vial contains peptide affinity-purified polyclonal antibody.
Synonyms:	L-PGDS, Lipocalin-PGDS
Immunogen:	Synthetic peptide from the internal region of human lipocalin-type PGDS
Cross Reactivity:	(-) H-PGDS
<b>Species Reactivity:</b>	(+) Human and mouse; other species not tested
Uniprot No.:	P41222
Form:	Lyophilized
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	TBS, pH 7.4, when reconstituted in 500 $\mu$ l deionized water
Host:	Rabbit
Application:	Western blot (WB); the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: Human CSF (20  $\mu$ g) Lane 2: Human CSF (10  $\mu$ g)

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, 11/14/2023

### 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

Prostaglandin D synthase (PGDS) catalyzes the isomerization of PGH<sub>2</sub> to produce PGD<sub>2</sub>. PGD<sub>2</sub> induces sleep, regulates nociception, inhibits platelet aggregation, and acts as an allergic mediator. Two distinct types of PGDS have been identified, namely the lipocalin-type enzyme ( $\beta$ -trace) and the hematopoietic enzyme.<sup>2,3</sup> Lipocalin-type PGDS is localized in the central nervous system and male genital organs of various mammals and the human heart whereas the hematopoietic PGDS is widely distributed in the peripheral tissues and localized in the antigen-presenting cells, mast cells, and megakaryocytes.<sup>2,4-6</sup> The lipocalin-type PGDS has been identified as  $\beta$ -trace, which is a major protein in human cerebrospinal fluid.<sup>3,7</sup> Human lipocalin-type PGDS is a 190 amino acid protein with a calculated molecular weight of 21,016, (observed at 24-26 kDa on immunoblot).<sup>1</sup>

#### References

- 1. Nagata, A., Suzuki, Y., Igarashi, M., et al. Proc. Natl. Acad. Sci. USA 88, 4020-4024 (1991).
- 2. Urade, Y., Watanabe, K., and Hayaishi, O. J. Lipid Mediators Cell Signalling 12, 257-273 (1995).
- 3. Toh, H., Kubodera, H., Nakajima, N., et al. Protein Engineering 9, 1067-1082 (1996).
- 4. Tokugawa, Y., Kunishige, I., Kuboto, Y., et al. Biol. Reprod. 58, 600-607 (1998).
- 5. Melegos, D.N., Diamandis, E.P., Oda, H., et al. Clin. Chem. 42, 1984-1991 (1996).
- 6. Kanaoka, Y., Fujimori, K., Kikuno, R., et al. Eur. J. Biochem. 267, 3315-3322 (2000).
- 7. Zahn, M., Mäder, A., Schmidt, B., et al. Neurosci. Lett. 154, 93-95 (1993).

#### 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