

PRODUCT INFORMATION



6-Thioguanosine-5'-O-diphosphate (sodium salt)

Item No. 38389

Formal Name: 6-thio-guanosine 5'-(trihydrogen diphosphate), trisodium salt

Synonyms: TGDP, 6-Thio-GDP

MF: C₁₀H₁₂N₅O₁₀P₂S • 3Na

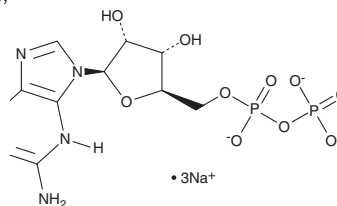
FW: 525.2

Purity: ≥95%

Supplied as: A solution in water

Storage: -80°C

Stability: ≥2 years



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

Description

6-Thioguanosine-5'-O-diphosphate is an inactive metabolite of the purine analog 6-thioguanine (Item No. 15774) and the immunosuppressive purine analog azathioprine (Item No. 13987).^{1,2} It is formed from 6-thioguanine *via* a 6-thioguanosine monophosphate intermediate by phosphorylation and from azathioprine *via* 6-mercaptopurine (Item No. 23675) and other intermediates by a non-enzymatic process and several enzymatic reactions. Elevated blood levels of 6-thioguanosine-5'-O-diphosphate in patients with Crohn's disease correlate with a lower response rate to azathioprine.²

References

1. Hawwa, A.F., Millership, J.S., Collier, P.A., *et al.* Pharmacogenomic studies of the anticancer and immunosuppressive thiopurines mercaptopurine and azathioprine. *Br. J. Clin. Pharmacol.* **66(4)**, 517-528 (2008).
2. Neurath, M.F., Kiesslich, R., Teichgraber, U., *et al.* 6-Thioguanosine diphosphate and triphosphate levels in red blood cells and response to azathioprine therapy in Crohn's disease. *Clin. Gastroenterol. Hepatol.* **3(10)**, 1007-1014 (2005).

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, 05/03/2023

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