

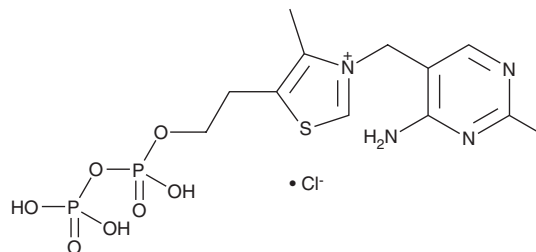
PRODUCT INFORMATION



Thiamine Pyrophosphate (chloride)

Item No. 20254

CAS Registry No.:	154-87-0
Formal Name:	3-[(4-amino-2-methyl-5-pyrimidinyl)methyl]-4-methyl-5-(4,6,6-trihydroxy-4,6-dioxido-3,5-dioxo-4,6-diphosphahex-1-yl)-thiazolium, monochloride
Synonyms:	Coccarboxylase, TPP, Vitamin B ₁ Pyrophosphate
MF:	C ₁₂ H ₁₉ N ₄ O ₇ P ₂ S • Cl
FW:	460.8
Purity:	≥95%
UV/Vis.:	λ _{max} : 246 nm
Supplied as:	A crystalline solid
Storage:	-20°C
Stability:	≥4 years



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

Laboratory Procedures

Thiamine pyrophosphate (TPP) (chloride) is supplied as a crystalline solid. Aqueous solutions of TPP (chloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of TPP (chloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

TPP (chloride) is a metabolite of vitamin B₁. It is a cofactor synthesized in the cytosol that is required for the activity of cytosolic transketolase and mitochondrial pyruvate, oxoglutarate, and branched-chain keto acid dehydrogenases.¹ TPP (chloride) prevents hyperglycemia-induced retinopathy and desflurane-induced hepatotoxicity in rats.^{2,3}

References

1. Hawkins, C.F., Borges, A., and Perham, R.N. A common structural motif in thiamin pyrophosphate-binding enzymes. *FEBS Lett.* **255(1)**, 77-82 (1989).
2. Cinici, E., Ahiskali, I., Cetin, N., *et al.* Effect of thiamine pyrophosphate on retinopathy induced by hyperglycemia in rats: A biochemical and pathological evaluation. *Indian J. Ophthalmol.* **64(6)**, 434-439
3. Arslan, A., Kuyrukluylidiz, U., Binici, O., *et al.* Can thiamine pyrophosphate prevent desflurane induced hepatotoxicity in rats? *Acta Cir. Bras.* **31(3)**, 168-175 (2016).

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, 12/21/2022

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