PRODUCT INFORMATION



Rp-8-pCPT-Cyclic GMPS (sodium salt)

Item No. 18445



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

Laboratory Procedures

Rp-8-pCPT-Cyclic GMPS (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the Rp-8-pCPT-cyclic GMPS (sodium salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Rp-8-pCPT-cGMPS is a stable, cell-permeable cGMP analog that competitively inhibits cGMP-dependent protein kinases (cGKs), including cGK I α and cGK II (IC₅₀s = 18.3 and 0.16 μ M, respectively).¹ The pCPT (p-chlorophenylthio) group at the 8-position of the purine increases both enzyme affinity and membrane permeability over related compounds.² At 10 μ M, it blocks the relaxation of rat tail arteries induced by the nitric oxide donor SIN-1 (Item No. 82220).³

References

- 1. Gamm, D.M., Francis, S.H., Angelotti, T.P., et al. The type II isoform of cGMP-dependent protein kinase is dimeric and possesses regulatory and catalytic properties distinct from the type I isoforms. J. Biol. Chem. 270(45), 27380-27388 (1995).
- 2. Kramer, R.H. and Tibbs, G.R. Antagonists of cyclic nucleotide-gated channels and molecular mapping of their site of action. J. Neurosci. 16(4), 1285-1293 (1996).
- Ouedraogo, S., Tschöp, M., Stoclet, J.-C., et al. Effects of cyclic GMP and analogues on neurogenic 3. transmission in the rat tail artery. Br. J. Pharmacol. 112, 867-872 (1994).

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

subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information luyer agrees to purchase the material can be found on our website

Copyright Cayman Chemical Company, 10/05/2022

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