

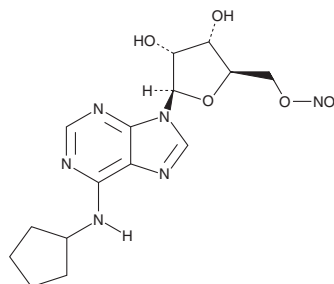
PRODUCT INFORMATION



INNO-8875

Item No. 33345

CAS Registry No.: 871108-05-3
Formal Name: N-cyclopentyl-adenosine, 5'-nitrate
Synonyms: PJ 875, Trabodenoson
MF: C₁₅H₂₀N₆O₆
FW: 380.4
Purity: ≥95%
UV/Vis.: λ_{max}: 268 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

INNO-8875 is supplied as a crystalline solid. A stock solution may be made by dissolving the INNO-8875 in the solvent of choice, which should be purged with an inert gas. INNO-8875 is soluble in the organic solvent DMSO.

Description

INNO-8875 is an adenosine A₁ receptor agonist.^{1,2} It binds to adenosine A₁ receptors (K_i = 0.97 nM) and is greater than 10,000-fold selective for adenosine A₁ over A₂ receptors. Topical application of INNO-8875 (25 µg/eye) reduces intraocular pressure in a rabbit model of ocular hypertension.² INNO-8875 (1-50 µg/kg) decreases atrial refractoriness and heart rate, as well as slows atrioventricular nodal conduction, in anesthetized rats.¹

References

1. Mor, M., Shalev, A., Dror, S., *et al.* INO-8875, a highly selective A₁ adenosine receptor agonist: Evaluation of chronotropic, dromotropic, and hemodynamic effects in rats. *J. Pharmacol. Exp. Ther.* **344**(1), 59-67 (2012).
2. Kim, N., Crosson, C., Supuran, C., *et al.* INO-8875, an adenosine A₁ agonist, in development for open-angle glaucoma reduces IOP in three rabbit models. *Invert. Ophthalmol. Vis. Sci.* **50**(13), 4061 (2009).

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, 11/03/2022

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