

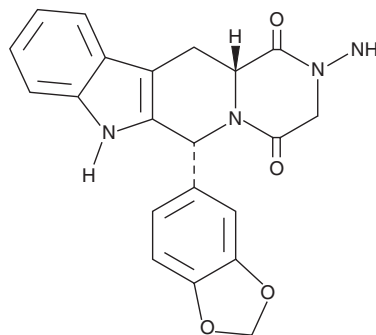
PRODUCT INFORMATION



Amino Tadalafil

Item No. 20876

CAS Registry No.: 385769-84-6
Formal Name: (6R,12aR)-2-amino-6-(1,3-benzodioxol-5-yl)-2,3,6,7,12,12a-hexahydro-pyrazino[1',2':1,6]pyrido[3,4-b]indole-1,4-dione
MF: C₂₁H₁₈N₄O₄
FW: 390.4
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 283, 290 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

Amino tadalafil is supplied as a crystalline solid. A stock solution may be made by dissolving the amino tadalafil in the solvent of choice, which should be purged with an inert gas. Amino tadalafil is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of amino tadalafil in these solvents is approximately 15 mg/ml.

Amino tadalafil is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, amino tadalafil should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Amino tadalafil has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Amino tadalafil is an analog of tadalafil (Item No. 14024), a potent inhibitor of phosphodiesterase 5 with applications in several conditions, including erectile dysfunction, pulmonary arterial hypertension, and lower urinary tract dysfunction.¹ Amino tadalafil has been detected as an illegal adulterant in a dietary supplement marketed for "tonic effect."²

References

1. Kim, D.-K., Ryu, D.H., Lee, N., *et al.* Synthesis and phosphodiesterase 5 inhibitory activity of new 5-phenyl-1,6-dihydro-7H-pyrazolo[4,3-d]pyrimidin-7-one derivatives containing an N-acylamido group phenyl ring. *Bioorg. Med. Chem.* **9(7)**, 1895-1899 (2001).
2. Hasegawa, T., Saijo, M., Ishi, T., *et al.* Structural elucidation of a tadalafil analogue found in a dietary supplement. *Shokuhin Eiseigaku Zasshi* **49(4)**, 311-315 (2008).

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, 11/10/2022

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