

PRODUCT INFORMATION

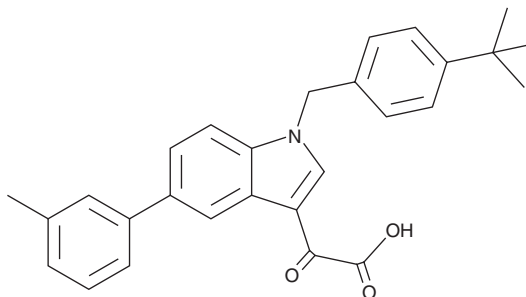


PAZ-417

Item No. 42122

CAS Registry No.: 481629-87-2
Formal Name: 1-[[4-(1,1-dimethylethyl)phenyl]methyl]-5-(3-methylphenyl)- α -oxo-1H-indole-3-acetic acid

Synonym: Aleplasinin
MF: C₂₈H₂₇NO₃
FW: 425.5
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

PAZ-417 is supplied as a solid. A stock solution may be made by dissolving the PAZ-417 in the solvent of choice, which should be purged with an inert gas. PAZ-417 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO

Description

PAZ-417 is an inhibitor of plasminogen activator inhibitor 1 (PAI-1; IC₅₀ = 655 nM).¹ It is selective for PAI-1 over tissue-type plasminogen activator (tPA) and plasminogen, as well as a panel of 40 neurotransmitter, endocrine, and enzymatic targets, at 10 μ M. PAZ-417 (5 μ M) enhances cleavage of monomeric and oligomeric amyloid- β (1-42) (A β 42) in a cell-free assay. *In vivo*, PAZ-417 (20 mg/kg) reduces plasma A β 40 and brain A β 40 and A β 42 levels in the Tg2576 transgenic mouse model of Alzheimer's disease. It also reverses cognitive deficits in the same model when administered at doses of 30 and 100 mg/kg.

Reference

1. Jacobsen, J.S., Comery, T.A., Martone, R.L., *et al.* Enhanced clearance of A β in brain by sustaining the plasmin proteolysis cascade. *Proc. Natl. Acad. Sci. USA* **105(25)**, 8754-8759 (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, 10/21/2024

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