

PRODUCT INFORMATION



MDK8582

Item No. 37307

CAS Registry No.: 185298-58-2

Formal Name: 4-(2-amino-2-oxoethoxy)-2-ethyl- α -oxo-1-(phenylmethyl)-1H-indole-3-acetamide

Synonyms: hnps-PLA Inhibitor, hnps-Phospholipase A Inhibitor, Human Nonpancreatic Secretory Phospholipase A Inhibitor, Human Nonpancreatic Secretory PLA Inhibitor

MF: $C_{21}H_{21}N_3O_4$

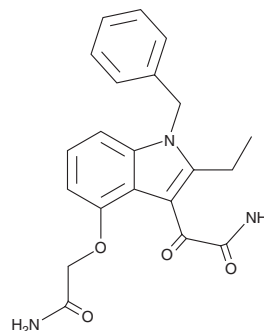
FW: 379.4

Purity: $\geq 98\%$

Supplied as: A 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

MDK8582 is supplied as a solid. A stock solution may be made by dissolving the MDK8582 in the solvent of choice, which should be purged with an inert gas. MDK8582 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

MDK8582 is an inhibitor of secretory phospholipase A_2 (sPLA $_2$; $IC_{50} = 0.124 \mu\text{M}$).¹

Reference

1. Bernard, P., Pintore, M., Berthon, J.Y., *et al.* A molecular modeling and 3D QSAR study of a large series of indole inhibitors of human non-pancreatic secretory phospholipase A_2 . *Eur. J. Med. Chem.* **36**(1), 1-19 (2001).

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, 04/26/2024

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