

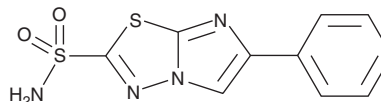
PRODUCT INFORMATION



AEG 3482

Item No. 26213

CAS Registry No.: 63735-71-7
Formal Name: 6-phenyl-imidazo[2,1-b]-1,3,4-thiadiazole-2-sulfonamide
MF: C₁₀H₈N₄O₂S₂
FW: 280.3
Purity: ≥98%
UV/Vis.: λ_{max}: 258 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

AEG 3482 is supplied as a crystalline solid. A stock solution may be made by dissolving the AEG 3482 in the solvent of choice. AEG 3482 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of AEG 3482 in ethanol is approximately 0.2 mg/ml and approximately 17 mg/ml in DMSO and DMF.

Description

AEG 3482 is an inhibitor of p75^{NTR}- and NRAGE-induced apoptosis.¹ It inhibits JNK signaling, caspase-3 cleavage, and cell death induced by the p75 neurotrophin receptor (p75^{NTR}) or the p75^{NTR}-interacting protein NRAGE in PC12^{rtTA} cells in a concentration-dependent manner. AEG 3482 inhibits cell death induced by withdrawal of nerve growth factor (NGF) in rat SCG neurons (EC₅₀ = ~20 μM). It also increases Hsp70 and Hsp25 levels in mouse embryonic fibroblasts (MEFs) in an HSF1-dependent manner and binds to recombinant Hsp90.

Reference

1. Salehi, A.H., Morris, S.J., Ho, W.C., *et al.* AEG3482 is an antiapoptotic compound that inhibits Jun kinase activity and cell death through induced expression of heat shock protein 70. *Chem. Biol.* **13(2)**, 213-223 (2006).

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, 12/20/2022

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