

PRODUCT INFORMATION



16F16

Item No. 17770

CAS Registry No.: 922507-80-0
Formal Name: 2-(2-chloroacetyl)-2,3,4,9-tetrahydro-1-methyl-1H-pyrido[3,4-b]indole-1-carboxylic acid, methyl ester

MF: C₁₆H₁₇ClN₂O₃

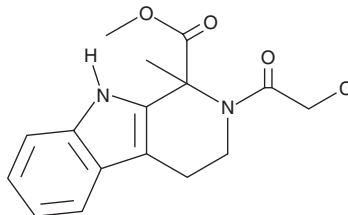
FW: 320.8

Purity: ≥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

16F16 is supplied as a solid. A stock solution may be made by dissolving the 16F16 in the solvent of choice, which should be purged with an inert gas. 16F16 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

16F16 is a protein disulfide isomerase (PDI) inhibitor.¹ It inhibits PDI reductase activity in an enzyme assay when used at concentrations ranging from 1 to 100 µg/ml.¹ 16F16 reduces PC12 cell apoptosis induced by the misfolded huntingtin protein HTT^{Q103}. It suppresses PDI-dependent mitochondrial outer membrane permeabilization (MOMP) in isolated PC12 cell mitochondria. 16F16 (2, 3, 4, and 10 µM) reduces HTT^{N90Q73} mutant huntingtin-induced medium spinal neuron death and MOMP in rat corticostriatal slices. It also reduces pyramidal neuron death induced by amyloid-β precursor protein (APP) in rat corticostriatal slices.

Reference

1. Hoffstrom, B.G., Kaplan, A., Letso, R., *et al.* Inhibitors of protein disulfide isomerase suppress apoptosis induced by misfolded proteins. *Nat. Chem. Biol.* **6(12)**, 900-906 (2010).

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, 12/19/2022

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