
PARP Inhibitor II, INH2BP

(Catalog No. M75003)

Alternative Names

5-Iodo-6-amino-1,2-benzopyrone

Description

A cell-permeable poly(ADP-ribose) polymerase (PARP) inhibitor. Offers protection against peroxynitrite and hydroxyl radicals in vitro and in vivo and abrogates peroxynitrite-induced mitochondrial transmembrane potential (DYm) reduction. Also exhibits malignant phenotype reversal in cancer cells and inhibits HIV-1 IIIb replication in vitro.

CAS Number

137881-27-7

Chemical Formula

C₉H₆INO₂

Molecular Weight

287.1

Size

5 mg

Purity

>98%

Solubility

DMSO

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
PARP Inhibitor II, INH2BP	5 mg	M75003 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.