

DPQ

(Catalog No. M71000)

Alternative Names

3,4-Dihydro-5-[4-(1-piperidinyl)butoxy]-1(2H)-isoquinolinone

Description

Cell-permeable. A very potent poly (ADP-ribose) polymerase (PARP) inhibitor

CAS Number

129075-73-6

Chemical Formula

$C_{18}H_{26}N_2O_2$

Molecular Weight

302.4

Size

1 mg

Purity

≥99% by NMR

Solubility

DMSO

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
DPQ	1 mg	M71000 -1

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