

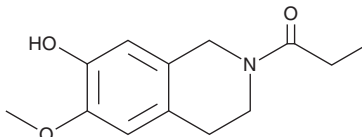
# PRODUCT INFORMATION



## PTIQ

Item No. 33540

**CAS Registry No.:** 1032822-42-6  
**Formal Name:** 1-(3,4-dihydro-7-hydroxy-6-methoxy-2(1H)-isoquinoliny)-1-propanone  
**MF:** C<sub>13</sub>H<sub>17</sub>NO<sub>3</sub>  
**FW:** 235.3  
**Purity:** ≥95%  
**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

PTIQ is supplied as a solid. A stock solution may be made by dissolving the PTIQ in the solvent of choice, which should be purged with an inert gas. PTIQ is slightly soluble in DMSO and methanol.

### Description

PTIQ is a neuroprotective compound.<sup>1</sup> It inhibits tetrahydro-L-biopterin-induced cytotoxicity and increases in matrix metalloproteinase-3 (MMP-3) levels in CATH.a cells (IC<sub>50</sub>s = 1.8 and 0.06 μM, respectively). PTIQ inhibits LPS-induced production of MMP-3, TNF-α, and COX-2 in BV-2 microglia (IC<sub>50</sub>s = 2.4, 6.5, and 9.3 μM, respectively). *In vivo*, PTIQ (3 and 30 mg/kg) increases survival of striatal nigral dopaminergic neurons, reduces microglial activation, and alleviates motor deficits in a mouse model of MPTP-induced Parkinson's disease.

### Reference

1. Son, H.J., Lee, J.A., Shin, N., *et al.* A novel compound PTIQ protects the nigral dopaminergic neurons in an animal model of Parkinson's disease induced by MPTP. *Br. J. Pharmacol.* **165**(7), 2213-2227 (2012).

#### 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, 11/15/2022

#### 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