

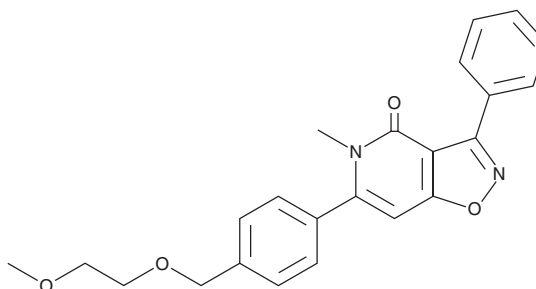
# PRODUCT INFORMATION



## IP7e

Item No. 29726

**CAS Registry No.:** 500164-74-9  
**Formal Name:** 6-[4-[(2-methoxyethoxy)methyl]phenyl]-5-methyl-3-phenylisoxazolo[4,5-c]pyridin-4(5H)-one  
**Synonym:** Isoxazolo-pyridinone 7e  
**MF:** C<sub>23</sub>H<sub>22</sub>N<sub>2</sub>O<sub>4</sub>  
**FW:** 390.4  
**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

IP7e is supplied as a solid. A stock solution may be made by dissolving the IP7e in the solvent of choice, which should be purged with an inert gas. IP7e is soluble in organic solvents such as ethanol (warmed) and DMSO. The solubility of IP7e in these solvents is approximately 20 and 100 mM, respectively.

### Description

IP7e is a brain-penetrant activator of nuclear receptor related protein 1 (Nurr1/NR4A2) signaling (EC<sub>50</sub> = 3.9 nM in a reporter assay).<sup>1</sup> *In vivo*, IP7e (10 mg/kg) prevents spinal cord axonal loss and demyelination and decreases spinal cord macrophage and T cell infiltration in a mouse model of experimental autoimmune encephalomyelitis (EAE).<sup>2</sup>

### References

1. Hintermann, S., Chiesi, M., von Krosigk, U., *et al.* Identification of a series of highly potent activators of the Nurr1 signaling pathway. *Bioorg. Med. Chem. Lett.* **17(1)**, 193-196 (2007).
2. Montarolo, F., Raffaele, C., Perga, S., *et al.* Effects of isoxazolo-pyridinone 7e, a potent activator of the Nurr1 signaling pathway, on experimental autoimmune encephalomyelitis in mice. *PLoS One* **9(9)**, e108791 (2014).

#### 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, 10/04/2022

#### 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