

SRPIN340

SRPK inhibitor

Catalog # S21-901

Lot # D3370-10

CAS # 218156-96-8

Product Description

Molecular Formula: C₁₈H₁₈F₃N₃O

Appearance: White powder

Molecular Weight: 349.4

Purity: >98% (TLC); NMR (Conforms)

Solubilization: May be dissolved in DMSO (40 mg/ml) or ethanol (25 mg/ml)

Alias

N-[2-(1-Piperidinyl)-5-(trifluoromethyl)phenyl]-4-pyridinecarboxamide

Specific Activity

SRPK inhibitor

Storage and Stability

Store desiccated as supplied at -20°C temperature for up to 2 years. Store solutions at -20°C for up to 1 month.

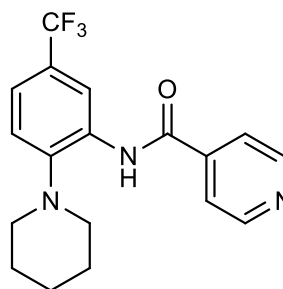
Scientific Background

Serine arginine protein kinase (SRPK) 1 inhibitor (K_i = 0.89 μM)¹. Inhibits SRPK2 at higher concentrations. Does not significantly inhibit other SRPKs such as CLK1 and CLK4, or other classes of SR kinases

References

1. T Fukuhara *et al. Proc. Natl. Acad. Sci. USA* 2006 103:11329
2. MV Gammons *et al. Invest. Ophthalmol. Vis. Sci.* 2013 54:5797
3. MV Gammons *et al. Br. J. Cancer* 2014 111:477

Molecular Structure



SRPIN340

SRPK inhibitor

Catalog Number	S21-901
CAS Number	218156-96-8
Specific Lot Number	D3370-10
Purity	>98% (TLC); NMR (Conforms)
Format	Lyophilized white powder
Solubilization	May be dissolved in DMSO (40 mg/ml) Ethanol (25 mg/ml)
Storage & Shipping	Store desiccated as supplied at -20°C temperature for up to 2 years. Store solutions at -20°C for up to 1 month.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.