

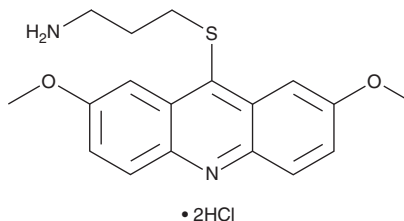
# PRODUCT INFORMATION



## LDN-192960 (hydrochloride)

Item No. 19089

**CAS Registry No.:** 2309172-48-1  
**Formal Name:** 3-[(2,7-dimethoxy-9-acridinyl)thio]-1-propanamine, dihydrochloride  
**MF:** C<sub>18</sub>H<sub>20</sub>N<sub>2</sub>O<sub>2</sub>S • 2HCl  
**FW:** 401.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 264 nm  
**Supplied as:** A crystalline 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

LDN-192960 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the LDN-192960 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. LDN-192960 (hydrochloride) is soluble in the organic solvent DMSO at a concentration of approximately 33 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of LDN-192960 (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of LDN-192960 (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

LDN-192960 is an inhibitor of haspin protein kinase and dual-specificity tyrosine-(Y)-phosphorylation regulated kinase 2 (DYRK2) with IC<sub>50</sub> values of 10 and 48 nM, respectively.<sup>1</sup> It is selective for haspin and DYRK2 over TRKB, CLK1, DYRK1A, DYRK3, ROS, HIPK1, HIPK2, PIM-1, and PIM-2 kinases (IC<sub>50</sub>'s = 720-91,000 nM).

### Reference

1. Cuny, G.D., Robin, M., Ulyanova, N.P., *et al.* Structure-activity relationship study of acridine analogs as haspin and DYRK2 kinase inhibitors. *Bioorg. Med. Chem. Lett.* **20(12)**, 3491-3493 (2010).

#### 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, 11/28/2022

#### 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