# **PRODUCT** INFORMATION



## HJC0152 (hydrochloride)

Item No. 22351

CAS Registry No.:	1420290-99-8	0 N
Formal Name:	2-(2-aminoethoxy)-5-chloro-N-(2-chloro-4-	U <sub>2</sub> N
	nitrophenyl)-benzamide, monohydrochloride	
MF:	$C_{15}H_{13}CI_2N_3O_4 \bullet HCI$	H • HCI
FW:	406.6	↓ N ↓ ↓
Purity:	≥98%	
UV/Vis.:	λ <sub>max</sub> : 326 nm	
Supplied as:	A crystalline solid	
Storage:	-20°C	H <sub>2</sub> N
Stability:	≥4 years	C C
Information represente	the product specifications. Batch specific analytical result	ts are provided on each certificate of analysis

#### Laboratory Procedures

HJC0152 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the HJC0152 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. HJC0152 (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of HJC0152 (hydrochloride) in these solvents is approximately 2 and 0.1 mg/ml, respectively.

#### Description

HJC0152 is an orally bioavailable inhibitor of STAT3 (32 and 62% inhibition at 10 and 20 µM, respectively, in MDA-MB-231 cells in a luciferase reporter assay).<sup>1</sup> It reduces total STAT3 and phosphorylated STAT3 levels in MDA-MB-231 cells and inhibits nuclear translocation of phosphorylated STAT3. HJC0152 inhibits proliferation of MCF-7 and MDA-MB-231 breast cancer and AsPC-1 and PANC-1 pancreatic cancer cells (IC<sub>50</sub>s = 0.91, 1.64, 1.9, and 1.08  $\mu$ M, respectively). It also halts the cell cycle at the G<sub>0</sub>/G<sub>1</sub> phase, induces apoptosis, and suppresses cell proliferation in human head and neck squamous cell carcinoma cells.<sup>2</sup> HJC0152 (7.5 mg/kg, i.p. or 25 mg/kg, p.o.) inhibits tumor growth in an MDA-MB-231 breast cancer mouse xenograft model and in an orthotopic mouse model of squamous cell carcinoma.<sup>1,2</sup>

#### References

- 1. Chen, H., Yang, Z., Ding, C., et al. Discovery of O-alkylamino tethered niclosamide derivatives as potent and orally bioavailable anticancer agents. ACS Med. Chem. Lett. 4(2), 180-185 (2013).
- 2. Wang, Y., Wang, S., Wu, Y., et al. Suppression of the growth and invasion of human head and neck squamous cell carcinomas via regulating STAT3 signaling and the miR-21/ $\beta$ -catenin axis with HJC0152. Mol. Cancer Ther. 16(4), 578-590 (2017).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFFTY 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

uyer 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, 11/14/2022

### 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