

# PRODUCT INFORMATION



## TP-238 (hydrochloride)

Item No. 27254

**CAS Registry No.:** 2415263-05-5  
**Formal Name:** 6-[4-[3-(dimethylamino)propoxy]phenyl]-2-(methylsulfonyl)-N-[3-(1H-pyrazol-1-yl)propyl]-4-pyrimidinamine, monohydrochloride

**MF:** C<sub>22</sub>H<sub>30</sub>N<sub>6</sub>O<sub>3</sub>S • HCl

**FW:** 495.0

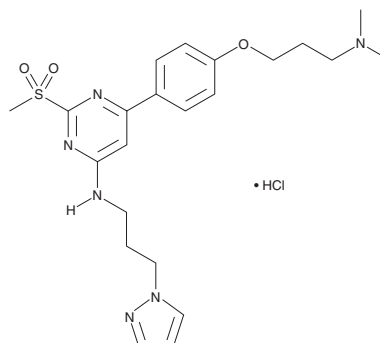
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 272 nm

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

TP-238 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the TP-238 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. TP-238 (hydrochloride) is soluble in the organic solvent DMSO at a concentration of approximately 1 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 TP-238 (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of TP-238 (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

TP-238 is a chemical probe for the bromodomains of cat eye syndrome chromosome region, candidate 2 (CECR2) and bromodomain PHD finger transcription factor (BPTF). It has on-target biochemical activity with CECR2 when used at concentrations ranging from 10 to 30 nM and with BPTF at concentrations ranging from 100 to 350 nM. It is selective for the CECR2 and BPTF bromodomains over BRD9 (IC<sub>50</sub> = 1.4 μM), as well as over a panel of 338 kinases at 1 μM. TP-238 binds to CECR2 and BPTF with K<sub>D</sub> values of 10 and 120 nM, respectively, in an isothermal titration calorimetry (ITC) assay and IC<sub>50</sub> values of 30 and 350 nM, respectively, in an AlphaScreen® assay. See the Structural Genomics Consortium (SGC) website for more information.

#### 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, 11/06/2020

#### 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