# **AZD1208**

#### pan-PIM kinase inhibitor

AZD1208 is a highly selective and potent pan-PIM kinase inhibitor. Displayed efficacy in models of acute myeloid leukemia, prostate cancer, and triple-negative breast cancer. Combination treatment of AZD1208 and imatinib eliminated chronic myelogenous leukemia stem cells in cell culture and a patient-derived mouse xenograft model. It improved immunotherapeutic antitumor T-cell response and disrupted the myeloid cell-mediated immunosuppressive tumor microenvironment increasing T-cell-mediated antitumor immunity.

## **Ordering Information**

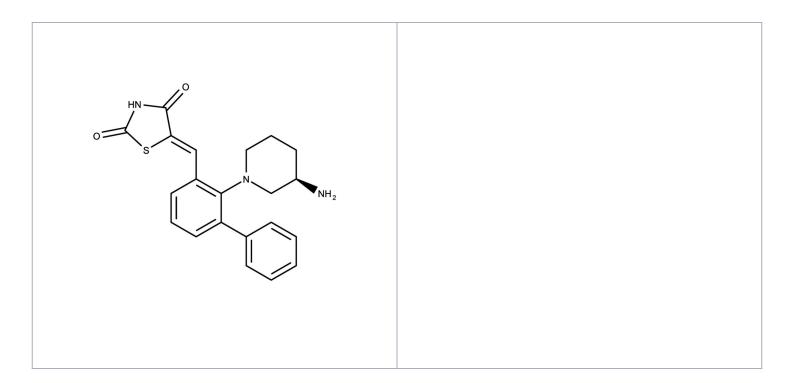
Order Online »

ENZ-CHM457-0025

25mg

Manuals, SDS & CofA

**View Online »** 



### **Handling & Storage**

**Use/Stability** As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

**Shipping** Ambient Temperature

# Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name (5Z)-5-[[2-[(3R)-3-Aminopiperidin-1-yl]-3-

phenylphenyl]methylidene]-1,3-thiazolidine-2,4-dione

Appearance Yellow solid.

CAS 1204144-28-4

Couple Type Inhibitor

 $\mathbf{C_{21}H_{21}N_{3}O_{2}S}$ 

**Identity** Determined by NMR.

MW 379.48

Purity ≥98% (HPLC)

**Solubility** Soluble in DMSO (2 mg/ml with warming).

Last modified: January 24, 2025