

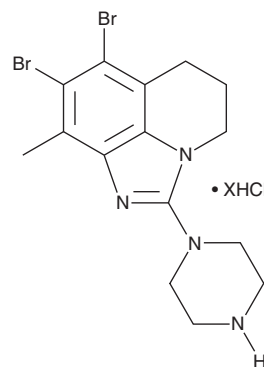
# PRODUCT INFORMATION



## SEL120-34A

Item No. 27906

**CAS Registry No.:** 1609452-30-3  
**Formal Name:** 7,8-dibromo-5,6-dihydro-9-methyl-2-(1-piperazinyl)-4H-imidazo[4,5,1-ij]quinoline, hydrochloride  
**MF:** C<sub>15</sub>H<sub>18</sub>Br<sub>2</sub>N<sub>4</sub> • XHCl  
**FW:** 414.1  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 228, 259, 300 nm  
**Supplied as:** A 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

SEL120-34A is supplied as a solid. A stock solution may be made by dissolving the SEL120-34A in the solvent of choice, which should be purged with an inert gas. SEL120-34A is soluble in organic solvents such as ethanol and DMSO.

### Description

SEL120-34A is a dual inhibitor of cyclin-dependent kinase 8 (Cdk8) and Cdk19 (IC<sub>50</sub>s = 4.4 and 10.4 nM, respectively).<sup>1</sup> It is selective for Cdk8 and Cdk19 over a panel of CDK/cyclin complexes at 1 μM. SEL120-34A (0.1-5 μM) inhibits IFN-γ-induced phosphorylation of STAT1 in HCT116 colon cancer cells and the proliferation of various acute myeloid leukemia (AML) cells, including MOLM-16, MV4-11, and KG-1 cells (GI<sub>50</sub>s = <1 μM). It reduces tumor growth in a KG-1 mouse xenograft model when administered at a dose of 60 mg/kg.

### Reference

1. Rzymiski, T., Mikula, M., Żytkiewicz, E., *et al.* SEL120-34A is a novel CDK8 inhibitor active in AML cells with high levels of serine phosphorylation of STAT1 and STAT5 transactivation domains. *Oncotarget* **8**(20), 33779-33795 (2017).

**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, 11/18/2022

### 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