
MLN8237 (Alisertib)

(Catalog No. M76000)

Alternative Names

Alisertib

Description

MLN8237 is a selective Aurora kinase A inhibitor with median IC₅₀ of 61 nM. MLN8237 (Alisertib) is a second-generation, orally bioavailable, highly selective small molecule inhibitor of the Aurora A kinase (serine/threonine protein kinase) with potential antineoplastic activity. MLN8237 (Alisertib) binds to and inhibits Aurora A kinase, which may result in disruption of the assembly of the mitotic spindle apparatus, disruption of chromosome segregation, and inhibition of cell proliferation.

CAS Number

1028486-01-2

Chemical Formula

C₂₇H₂₀ClFN₄O₄

Molecular Weight

518.92

Size

5 mg

Purity

99.27%

Solubility

DMSO ≥44mg/mL Water <1mg/mL Ethanol <1mg/mL

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
MLN8237 (Alisertib)	5 mg	M76000 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.