

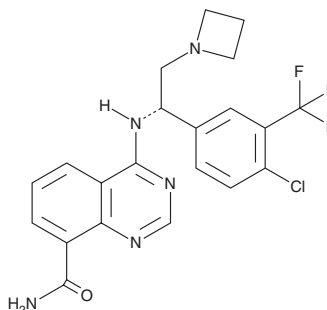
# PRODUCT INFORMATION



**M2698**

Item No. 36816

**CAS Registry No.:** 1379545-95-5  
**Formal Name:** 4-[[[(1S)-2-(1-azetidiny)-1-[4-chloro-3-(trifluoromethyl)phenyl]ethyl]amino]-8-quinazolinecarboxamide  
**Synonym:** MSC2363318A  
**MF:** C<sub>21</sub>H<sub>19</sub>ClF<sub>3</sub>N<sub>5</sub>O  
**FW:** 449.9  
**Purity:** ≥98%  
**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

M2698 is supplied as a solid. A stock solution may be made by dissolving the M2698 in the solvent of choice, which should be purged with an inert gas. M2698 is soluble in DMSO.

## Description

M2698 is a dual inhibitor of p70 ribosomal S6 kinase (p70S6K) and Akt1 (IC<sub>50</sub>s = 1.1 and 4 nM, respectively).<sup>1</sup> It is selective for p70S6K and Akt1 over Aurora B kinase (IC<sub>50</sub> = 170 nM), as well as a panel of 265 kinases at 1 μM, but also inhibits protein kinase A (PKA), p90 ribosomal S6 kinase 1 (MSK1), MSK2, cGMP-dependent protein kinase 1α (PKG1α), PKG1β, and protein kinase X (PRKX). M2698 reduces tumor volume in MiaPaCa-2 pancreatic cancer and A549 non-small cell lung cancer (NSCLC) mouse xenograft models when administered at doses of 20 or 30 mg/kg, respectively.

## Reference

1. DeSelm, L., Huck, B., Lan, R., *et al.* Identification of clinical candidate M2698, a dual p70S6K and Akt inhibitor, for treatment of PAM pathway-altered cancers. *J. Med. Chem.* **64**(19), 14603-14619 (2021).

### 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, 12/09/2022

## 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