

CERTIFICATE of ANALYSIS



FIIN-3

N-[4-[[[(2,6-dichloro-3,5-dimethoxyphenyl)amino]carbonyl][6-[[4-(4-methyl-1-piperazinyl)phenyl]amino]-4-pyrimidinyl]amino]methyl]phenyl]-2-propenamide

Item No. 20709 * Batch No. 0492667

Purity Specification: $\geq 98\%$

Molecular formula: C₃₄H₃₆Cl₂N₈O₄

CAS number: 1637735-84-2

Formula weight: 691.60

Expiration date: 19OCT2028

Overview

Tests

Results

HPLC	Purity: 98.0 %
IR	Conforms
Mass Spec	MH+: 691.4
Melting point	183 - 185 °C
TLC	Purity: 100 %
UV	λ max: 287.00 nm

Reviewed and approved by: Jennifer LaBrecque

WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

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, 10/12/2018

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