

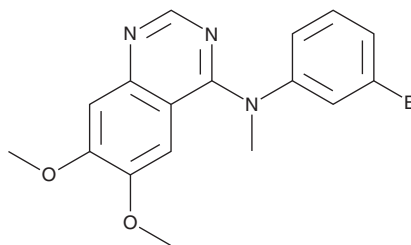
PRODUCT INFORMATION



EBE-A22

Item No. 43954

CAS Registry No.: 229476-53-3
Formal Name: N-(3-bromophenyl)-6,7-dimethoxy-N-methyl-4-quinazolinamine
MF: $C_{17}H_{16}BrN_3O_2$
FW: 374.2
Purity: $\geq 98\%$
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

EBE-A22 is supplied as a solid. A stock solution may be made by dissolving the EBE-A22 in the solvent of choice, which should be purged with an inert gas. EBE-A22 is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

EBE-A22 is an inactive derivative of PD 153035 (Item Nos. 18080 | 14879).¹

Reference

1. Bridges, A.J., Zhou, H., Cody, D.R., *et al.* Tyrosine kinase inhibitors. 8. An unusually steep structure-activity relationship for analogues of 4-(3-bromoanilino)-6,7-dimethoxyquinazoline (PD 153035), a potent inhibitor of the epidermal growth factor receptor. *J. Med. Chem.* **39**(1), 267-276 (1996).

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, 08/15/2025

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