

PRODUCT INFORMATION



UMB-32

Item No. 17123

CAS Registry No.: 1635437-39-6

Formal Name: N-(1,1-dimethylethyl)-2-[4-(3,5-dimethyl-4-isoxazolyl)phenyl]-imidazo[1,2-a]pyrazin-3-amine

MF: $C_{21}H_{23}N_5O$

FW: 361.4

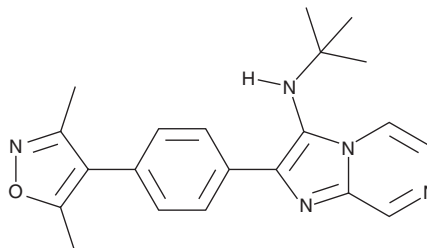
Purity: $\geq 98\%$

UV/Vis.: λ_{\max} : 258, 345 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Laboratory Procedures

UMB-32 is supplied as a crystalline solid. A stock solution may be made by dissolving the UMB-32 in the solvent of choice. UMB-32 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of UMB-32 in ethanol and DMF is approximately 20 mg/ml and approximately 10 mg/ml in DMSO.

UMB-32 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, UMB-32 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. UMB-32 has a solubility of approximately 0.1 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

UMB-32 is an inhibitor of the BET bromodomain BRD4 ($K_d = 550$ nM; $IC_{50} = 637$ nM) and the bromodomain-containing transcription factor TAF1 ($K_d = 560$ nM) and TAF1L ($K_d = 1.3$ μM).¹

Reference

1. McKeown, M.R., Shaw, D.L., Fu, H., *et al.* Biased multicomponent reactions to develop novel bromodomain inhibitors. *J. Med. Chem.* **57**(21), 9019-9027 (2014).

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, 07/21/2016

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