

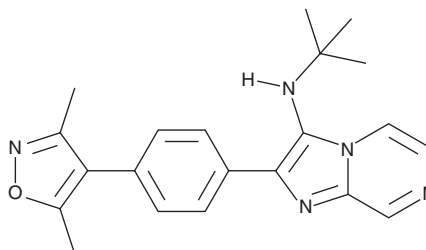
PRODUCT INFORMATION



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: C₂₁H₂₃N₅O
FW: 361.4
Purity: ≥98%
UV/Vis.: λ_{max}: 258, 345 nm
Supplied as: A crystalline 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

UMB-32 is supplied as a crystalline solid. A stock solution may be made by dissolving the UMB-32 in the solvent of choice, which should be purged with an inert gas. UMB-32 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). 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, 02/26/2025

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