

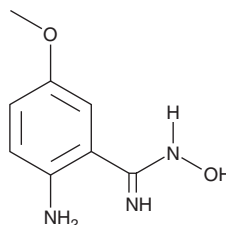
PRODUCT INFORMATION



para-methoxy-2-amino Benzamidoxime

Item No. 9004402

CAS Registry No.: 1359828-78-6
Formal Name: 2-amino-N-hydroxy-5-methoxy-benzenecarboximidamide
Synonyms: PMA, *p*-methoxy-2-amino Benzamidoxime
MF: C₈H₁₁N₃O₂
FW: 181.2
Purity: ≥95%
Ex./Em. Max: 389/515 nm
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

para-methoxy-2-amino-Benzamidoxime (PMA) is supplied as a solid. A stock solution may be made by dissolving the PMA in the solvent of choice, which should be purged with an inert gas. PMA is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in methanol and acetonitrile.

Description

PMA is a fluorescent and reactive probe for ketones.¹ Upon binding to ketones, PMA displays excitation/emission maxima of 389 and 515 nm, respectively. It has been used in the detection of alcohol dehydrogenases (ADHs), which convert alcohols to ketones.

Reference

1. Mei, Z., Zhang, K., Qu, G., *et al.* High-throughput fluorescence assay for ketone detection and its applications in enzyme mining and protein engineering. *ACS Omega* **5**(23), 13588-13594 (2020).

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, 05/07/2024

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