

PRODUCT INFORMATION



Methyl diaminobenzene-BODIPY (trifluoroacetate salt)

Item No. 44873

Formal Name: (T-4)-5-dimethyl-2H-pyrrol-2-ylidene-κN¹] methyl]-2,4-dimethyl-1Hpyrrole-3-carboxylato-κN] difluoro-[ethyl 5-[(4,5-diamino-2-methylphenyl) [4-(ethoxycarbonyl)-3 boron, 2,2,2-trifluoroacetate

Synonyms: MBo, Methylglyoxal Probe, Methylglyoxal Sensor, MGO Probe, MGO Sensor

MF: C₂₆H₃₁BF₂N₄O₄ • CF₃COOH

FW: 626.4

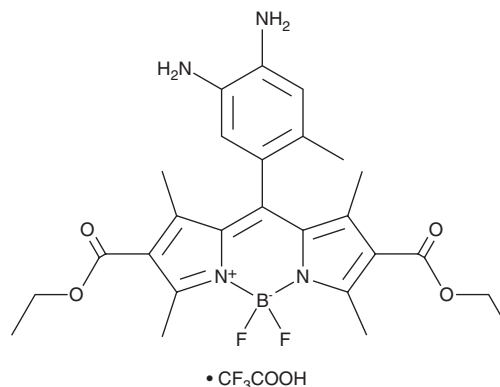
Purity: ≥95%

Abs./Em. Max: 496/530 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

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

Description

Methyl diaminobenzene-BODIPY is a turn-on fluorescent probe for the detection of methylglyoxal.¹ It displays absorbance/emission maxima of 496/530 nm, respectively, in DMSO and its fluorescence intensity increases in the presence of methylglyoxal. Methyl diaminobenzene-BODIPY is selective for methylglyoxal over glyoxal, glucosone, pyruvate, 1-phenyl-1,2-propanedione, phenylglyoxal, nitric oxide (NO) donors, and hydrogen peroxide. It has been used to quantify methylglyoxal levels in living cells and plasma.

Reference

1. Wang, T., Douglass, E.F., Jr., Fitzgerald, K.J., et al. A "turn-on" fluorescent sensor for methylglyoxal. *J. Am. Chem. Soc.* **135**(33), 12429-12433 (2013).

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, 01/20/2026

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