

PRODUCT INFORMATION

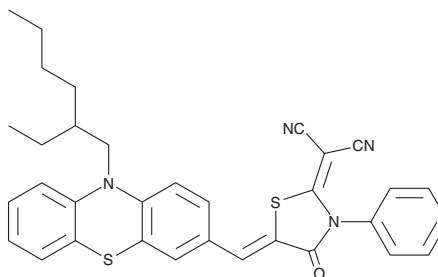


MIT-PZR

Item No. 28925

CAS Registry No.: 2376848-06-3
Formal Name: 2-[5-[[10-(2-ethylhexyl)-10H-phenothiazin-3-yl]methylene]-4-oxo-3-phenyl-2-thiazolidinylidene]-propanedinitrile

MF: C₃₃H₃₀N₄OS₂
FW: 562.7
Purity: ≥98%
UV/Vis.: λ_{max}: 242, 304, 352, 472 nm
Abs./Em. Max: 485/705 nm, respectively
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

MIT-PZR is supplied as a crystalline solid. A stock solution may be made by dissolving the MIT-PZR in the solvent of choice, which should be purged with an inert gas. MIT-PZR is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of MIT-PZR in these solvents is approximately 30 mg/ml.

Description

MIT-PZR is a mitochondria-targeted fluorescent probe.¹ It is aggregation-induced emission (AIE) active and displays absorption/emission maxima of 485/705 nm, respectively. MIT-PZR can be used in live cells and *in vivo*.

Reference

1. Dong, Y., Chen, Z., Hou, L., *et al.* Mitochondria-targeted aggregation-induced emission active near infrared fluorescent probe for real-time imaging. *Spectrochim. Acta. A. Mol. Biomol. Spectrosc.* **224**, 117456 (2019).

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, 05/21/2020

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