

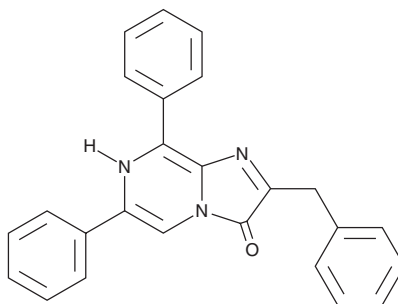
PRODUCT INFORMATION



Diphenylterazine

Item No. 37334

CAS Registry No.: 344940-63-2
Formal Name: 6,8-diphenyl-2-(phenylmethyl)-imidazo[1,2-a]pyrazin-3(7H)-one
Synonym: DTZ
MF: C₂₅H₁₉N₃O
FW: 377.4
Purity: ≥98%
UV/Vis.: λ_{max}: 259, 450 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

Diphenylterazine is supplied as a solid. A stock solution may be made by dissolving the diphenylterazine in the solvent of choice, which should be purged with an inert gas. Diphenylterazine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of diphenylterazine in DMSO and DMF is approximately 2 and 11 mg/ml, respectively. Diphenylterazine is slightly soluble in ethanol.

Description

Diphenylterazine is a synthetic bioluminescent luciferin that displays an emission maximum of 502 nm.¹ It has been used to quantify extracellular hydrogen peroxide when coupled to a boronated ester cage.²

References

1. Yeh, H.-W., Karmach, O., Ji, A., *et al.* Red-shifted luciferase-luciferin pairs for enhanced bioluminescence imaging. *Nat. Methods* **14**(10), 971-974 (2017).
2. O'Sullivan, J.J. and Heffrin, M.C. Development of an ATP-independent bioluminescent probe for detection of extracellular hydrogen peroxide. *Org. Biomol. Chem.* **20**(31), 6231-6238 (2022).

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, 03/07/2023

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