

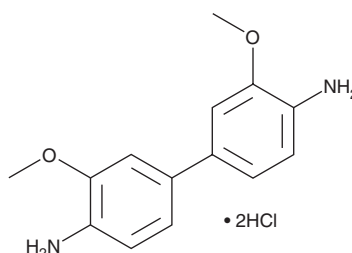
PRODUCT INFORMATION



o-Dianisidine (hydrochloride)

Item No. 17316

CAS Registry No.: 20325-40-0
Formal Name: 3,3'-dimethoxy-[1,1'-biphenyl]-4,4'-diamine, dihydrochloride
MF: C₁₄H₁₆N₂O₂ • 2HCl
FW: 317.2
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 297 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

o-Dianisidine (hydrochloride) is supplied as a crystalline solid. Aqueous solutions of *o*-dianisidine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of *o*-dianisidine (hydrochloride) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

o-Dianisidine (dihydrochloride) is a colorimetric peroxidase substrate that has been used as a substrate to semi-quantitatively measure lactate, uric acid (Item No. 16219), and glucose.^{1,2} It has also been used as a reagent for the detection of various metals, thiocyanates, and nitrites.³ This aromatic amine is initially colorless but turns to violet upon oxidization and is spectrophotometrically measured at a wavelength of 405 nm.

References

1. Sugiura, M. and Hirano, K. A new colorimetric method for determination of serum glucose. *Clin. Chim. Acta* **75(3)**, 387-391 (1977).
2. Mesulam, M.M. and Rosene, D.L. Sensitivity in horseradish peroxidase neurohistochemistry: A comparative and quantitative study of nine methods. *J. Histochem. Cytochem.* **27(3)**, 763-773 (1979).
3. National Toxicology Program. 3,3'-Dimethoxybenzidine and dyes metabolized to 3,3'-dimethoxybenzidine. *Rep. Carcinog.* **12**, 164-165 (2011).

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, 10/25/2022

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