

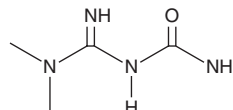
PRODUCT INFORMATION



Metformin hydroxy analog 2

Item No. 36316

CAS Registry No.: 98026-16-5
Formal Name: N-[(dimethylamino)iminomethyl]-urea
Synonym: N-Carbonylguanidine
MF: C₄H₁₀N₄O
FW: 130.1
Purity: ≥90%
UV/Vis.: λ_{max}: 229 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

Metformin hydroxy analog 2 is supplied as a solid. A stock solution may be made by dissolving the metformin hydroxy analog 2 in the solvent of choice, which should be purged with an inert gas. Metformin hydroxy analog 2 is soluble in the organic solvent dimethyl formamide at a concentration of approximately 10 mg/ml. Metformin hydroxy analog 2 is slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of metformin hydroxy analog 2 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of metformin hydroxy analog 2 in PBS (pH 7.2) is approximately 0.09 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Metformin hydroxy analog 2 is an oxidation product of the biguanide metformin (Item No. 13118).¹

Reference

1. Gholivand, M.B. and Mohammadi-Behzad, L. Differential pulse voltammetric determination of metformin using copper-loaded activated charcoal modified electrode. *Anal. Biochem.* **38(1)**, 53-60 (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, 11/07/2022

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