

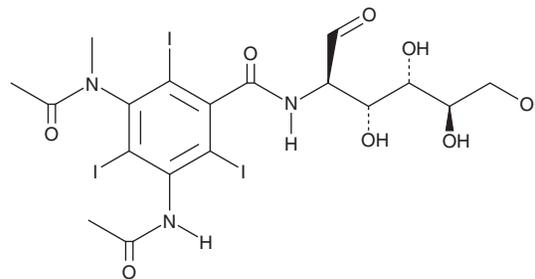
PRODUCT INFORMATION



Metrizamide

Item No. 27724

CAS Registry No.: 31112-62-6
Formal Name: 2-[[3-(acetylamino)-5-(acetylmethylamino)-2,4,6-triiodobenzoyl]amino]-2-deoxy-D-glucose
MF: C₁₈H₂₂I₃N₃O₈
FW: 789.1
Purity: ≥95%
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

Metrizamide is supplied as a solid. A stock solution may be made by dissolving the metrizamide in the solvent of choice, which should be purged with an inert gas. Metrizamide is soluble in organic solvents such as methanol and DMSO.

Description

Metrizamide is a non-ionic contrast agent.¹ *In vivo*, metrizamide (0.1 ml/kg of a 370 mg l/ml solution) does not induce EEG abnormalities, indicating minimal neurotoxicity in rabbits. It has also been used as a solute in gradient centrifugation applications.²

References

1. Maly, P., Elmqvist, D., Almén, T., *et al.* Comparison between EEG and observation of rabbit behaviour in evaluation of subarachnoid neurotoxicity of metrizamide. *Acta. Radiol. Diagn. (Stockh.)* **27(2)**, 235-240 (1986).
2. Hsu, C.-H., Di Carlo, D., Chen, C., *et al.* Microvortex for focusing, guiding and sorting of particles. *Lab Chip.* **8(12)**, 2128-2134 (2008).

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, 12/08/2022

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