

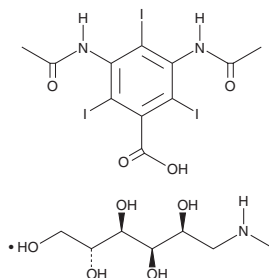
PRODUCT INFORMATION



Diatrizoate (meglumine)

Item No. 44267

CAS Registry No.: 131-49-7
Formal Name: 1-deoxy-1-(methylamino)-D-glucitol, 3,5-bis(acetylamino)-2,4,6-triiodobenzoate (1:1)
MF: $C_{11}H_9I_3N_2O_4 \bullet C_7H_{17}NO_5$
FW: 809.1
Purity: $\geq 98\%$
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

Diatrizoate (meglumine) is supplied as a solid. Aqueous solutions of diatrizoate (meglumine) can be prepared by directly dissolving the solid in aqueous buffers. Diatrizoate (meglumine) is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Diatrizoate is an iodinated and radiopaque contrast agent.^{1,2} Formulations containing diatrizoate have been used as radiographic contrast agents in gastrointestinal or urinary tract examinations.

References

1. Almén, T. and Trägårdh, B. Effects of non-ionic contrast media on the blood flow through the femoral artery of the dog. *Acta Radiol. Suppl.* **335**, 197-202 (1973).
2. Golman, K. and Almén, T. Contrast media-induced nephrotoxicity. Survey and present state. *Invest. Radiol.* **20**(1 Suppl.), S92-S97 (1985).

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, 10/06/2025

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