

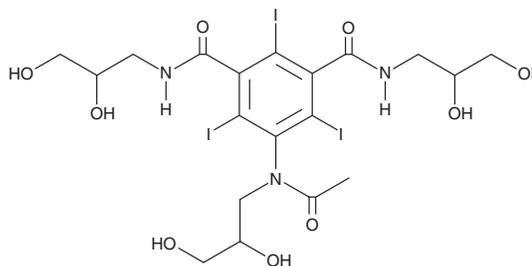
PRODUCT INFORMATION



Iohexol

Item No. 23753

CAS Registry No.: 66108-95-0
Formal Name: 5-[acetyl(2,3-dihydroxypropyl)amino]-N¹,N³-bis(2,3-dihydroxypropyl)-2,4,6-triiodo-1,3-benzenedicarboxamide
MF: C₁₉H₂₆I₃N₃O₉
FW: 821.1
Purity: ≥98% (mixture of isomers)
UV/Vis.: λ_{max}: 245 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

Iohexol is supplied as a crystalline solid. A stock solution may be made by dissolving the iohexol in the solvent of choice, which should be purged with an inert gas. Iohexol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of iohexol in ethanol is approximately 2 mg/ml and approximately 10 mg/ml in DMSO and DMF.

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 iohexol can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of iohexol in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Iohexol is a non-ionic and water-soluble contrast agent.¹ It induces inhibition of electrical activity but has no excitatory effect *in vitro* in rat hippocampal slices and lacks any effect *in vivo* on ventral root flexes of feline spinal cord, indicating minimal neurotoxicity. Iohexol also induces a smaller decrease in contractile force than the ionic and high osmolarity contrast agent metrizoate in isolated rabbit hearts perfused with a 350 mg/ml solution.² Formulations containing iohexol have been used for visualization of arteries, veins, the urinary tract, and joints using X-ray and computed tomography (CT) scan imaging techniques.

References

1. Bryan, R.N., Centeno, R.S., Hershkowitz, N., *et al.* Neurotoxicity of iohexol: A new nonionic contrast medium. *Radiology* **145**(2), 379-382 (1982).
2. Almén, T. and Bååth, L. Effects of iopentol, iohexol and metrizoate on the contractility of the isolated rabbit heart. *Acta Radiol. Suppl.* **370**, 61-63 (1987).

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, 11/29/2022

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