

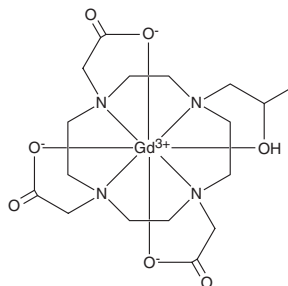
PRODUCT INFORMATION



Gadoteridol

Item No. 33577

CAS Registry No.: 120066-54-8
Formal Name: [10-[2-(hydroxy-κO)propyl]-1,4,7,10-tetraazacyclododecane-1,4,7-triacetato(3-)-κN¹,κN⁴,κN⁷,κN¹⁰,κO¹,κO⁴,κO⁷]-gadolinium
Synonyms: Gd-HP-D03A, SQ 32,692
MF: C₁₇H₂₉GdN₄O₇
FW: 558.7
Purity: ≥90%
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

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

Description

Gadoteridol is a gadolinium-based MRI contrast agent.^{1,2} It enhances the visibility of intraparenchymal and meningeal tumors in an orthotopic rat model of brain cancer when administered at a dose of 0.25 mmol/kg.¹ *In vitro* labeling of murine mesenchymal stem cells (MSCs) with gadoteridol and subsequent injection into mice has been used to track their migration to spinal cord injury lesions *in vivo*.² Formulations containing gadoteridol have been used to visualize intra- and extracranial lesions in the head, neck, and spine.

References

1. Runge, V.M., Gelblum, D.Y., and Jacobson, S. Gd HP-DO3A—experimental evaluation in brain and renal MR. *Magn. Reson. Imaging* **9**(1), 79-87 (1991).
2. Filippi, M., Boido, M., Pasquino, C., *et al.* Successful *in vivo* MRI tracking of MSCs labeled with gadoteridol in a spinal cord injury experimental model. *Exp. Neurol.* **282**, 66-77 (2016).

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, 11/17/2022

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