

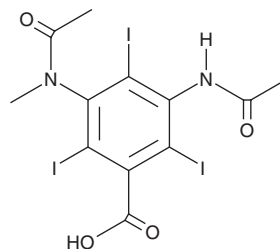
# PRODUCT INFORMATION



## Metrizoic acid

Item No. 38512

CAS Registry No.: 1949-45-7  
Formal Name: 3-(acetilamino)-5-(acetilmethylamino)-  
2,4,6-triiodo-benzoic acid  
MF:  $C_{12}H_{11}I_3N_2O_4$   
FW: 627.9  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^\circ\text{C}$   
Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Metrizoic acid is supplied as a solid. A stock solution may be made by dissolving the metrizoic acid in the solvent of choice, which should be purged with an inert gas. Metrizoic acid is soluble in methanol.

### Description

Metrizoic acid is an iodinated monocarboxylic acid.<sup>1</sup> It has been used in gradient media to isolate the parasite *B. hominis* from human feces as well as in live lymphocytes in adoptive transfer experiments in mice.<sup>1,2</sup> A salt form of metrizoic acid, metrizoate, previously has been used as an ionic contrast agent for X-ray imaging and angiography.<sup>3</sup>

### References

1. Lanuza, M.D., Carbajal, J.A., Villar, J., et al. Description of an improved method for *Blastocystis hominis* culture and axenization. *Parasitol. Res.* **83**(1), 60-63 (1996).
2. Korbely, M. and Dougherty, G.J. Photodynamic therapy-mediated immune response against subcutaneous mouse tumors. *Cancer Res.* **59**(8), 1941-1946 (1999).
3. 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, 01/19/2024

#### 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