

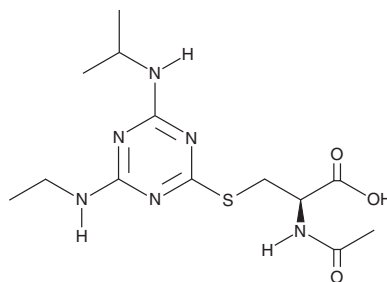
PRODUCT INFORMATION



Atrazine Mercapturate

Item No. 20395

CAS Registry No.: 138722-96-0
Formal Name: N-acetyl-S-[4-(ethylamino)-6-[(1-methylethyl)amino]-1,3,5-triazin-2-yl]-L-cysteine
Synonym: ATR Mercapturate
MF: C₁₃H₂₂N₆O₃S
FW: 342.4
Purity: ≥98%
UV/Vis.: λ_{max}: 223 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

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

Description

Atrazine mercapturate is a metabolite of the herbicide atrazine (Item No. 13375).¹ Atrazine has been reported to cause cancer in certain laboratory animals, have diverse effects on the reproductive system in animals and humans, and disrupt the normal function of the endocrine system.²⁻⁴ Atrazine mercapturate is readily measured in urine samples.^{1,5}

References

1. Koivunen, M.E., Dettmer, K., Vermeulen, R., *et al.* Improved methods for urinary atrazine mercapturate analysis—assessment of an enzyme-linked immunosorbent assay (ELISA) and a novel liquid chromatography-mass spectrometry (LC-MS) method utilizing online solid phase extraction (SPE). *Anal. Chim. Acta* **572(2)**, 180-189 (2006).
2. Eldridge, J.C., Wetzel, L.T., Stevens, J.T., *et al.* The mammary tumor response in triazine-treated female rats: A threshold-mediated interaction with strain and species-specific reproductive senescence. *Steroids* **64**, 672-678 (1999).
3. Goldman, J.M., Stoker, T.E., and Cooper, R.L. Neuroendocrine and reproductive effects of contemporary-use pesticides. *Toxicol. Ind. Health* **15**, 26-36 (1999).
4. Hayes, T.B., Collins, A., Lee, M., *et al.* Hermaphroditic, demasculinized frogs after exposure to the herbicide atrazine at low ecologically relevant doses. *Proc. Natl. Acad. Sci. USA* **99(8)**, 5476-5480 (2002).
5. Chevrier, C., Limon, G., Monfort, C., *et al.* Urinary biomarkers of prenatal atrazine exposure and adverse birth outcomes in the PELAGIE birth cohort. *Environ. Health Perspect.* **119(7)**, 1034-1041 (2011).

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, 09/11/2023

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