

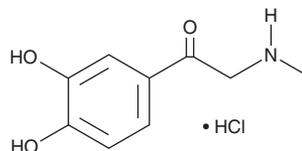
PRODUCT INFORMATION



Adrenalone (hydrochloride)

Item No. 43563

CAS Registry No.: 62-13-5
Formal Name: 1-(3,4-dihydroxyphenyl)-2-(methylamino)-ethanone, monohydrochloride
MF: C₉H₁₁NO₃ • HCl
FW: 217.7
Purity: ≥95%
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

Adrenalone (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the adrenalone (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Adrenalone (hydrochloride) is soluble (≥10 mg/ml) in DMSO.

Description

Adrenalone is an oxidized derivative of epinephrine.¹ It induces dsDNA breaks in RFL-6 rat fetal lung fibroblasts when used at a concentration of 250 μM. Adrenalone (20 mg/kg) increases the blood levels of fibrinogen in rats.² Formulations containing adrenalone have previously been used as topical agents to prolong the activity of anesthetics.

References

1. Murakami, H., Yamada, K., Shirahata, S., *et al.* Double strand breakage of chromatin DNA in cultured mammalian cells and DNA extracted from λ phage by aromatic reductones derived from adrenalone. *Agric. Biol. Chem.* **42(5)**, 1019-1023 (1978).
2. Henriques, S.B., Henriques, O.B., and Levy, A. Slow hyperfibrinogenaemic action of adrenaline and related substances. *Br. J. Pharmacol. Chemother.* **11(2)**, 99-103 (1956).

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, 06/20/2025

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