

PRODUCT INFORMATION



Acid Violet 17

Item No. 44936

CAS Registry No.: 4129-84-4
Formal Name: N-[4-[[4-(diethylamino)phenyl][4-[ethyl[(3-sulphophenyl)methyl]amino]phenyl]methylene]-2,5-cyclohexadien-1-ylidene]-N-ethyl-3-sulfo-benzenemethanaminium, inner salt, monosodium salt

Synonyms: AV17, C.I. 42640, Coomassie Violet R200

MF: $C_{41}H_{44}N_3O_6S_2 \cdot Na$

FW: 761.9

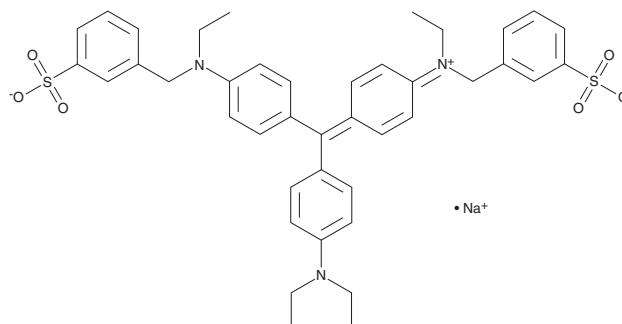
Absorbance: 548 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Special Conditions: Light sensitive



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

Laboratory Procedures

Acid violet 17 is supplied as a solid. A stock solution may be made by dissolving the acid violet 17 in the solvent of choice, which should be purged with an inert gas. Acid violet 17 is soluble (≥ 10 mg/ml) in DMSO.

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 acid violet 17 can be prepared by directly dissolving the solid in aqueous buffers. Acid violet 17 is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Acid violet 17 is an anionic dye.¹ It displays an absorbance maximum at 548 nm. Acid violet 17 has been found as a contaminant in textile wastewaters. Formulations containing acid violet 17 have been used as textile dyes.

Reference

1. Sharma, D.K., Saini, H.S., Singh, M., *et al.* Biological treatment of textile dye acid violet-17 by bacterial consortium in an up-flow immobilized cell bioreactor. *Let. Appl. Microbiol.* **38**(5), 345-350 (2004).

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, 02/06/2026

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