

# PRODUCT INFORMATION

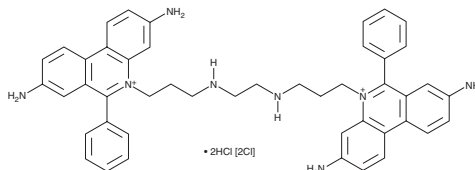


## Ethidium Homodimer-1

Item No. 43850

**CAS Registry No.:** 61926-22-5  
**Formal Name:** 5,5'-[1,2-ethanediylbis(imino-3,1-propanediyl)]bis[3,8-diamino-6-phenyl-phenanthridinium, dihydrochloride, dichloride

**Synonym:** EthD-1  
**MF:** C<sub>46</sub>H<sub>48</sub>N<sub>8</sub> • 2HCl [2Cl]  
**FW:** 856.8  
**Abs./Em. Max:** 528/617 nm  
**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.

### Description

Ethidium homodimer-1 is a membrane-impermeable fluorescent dye and DNA intercalator.<sup>1,2</sup> It displays absorbance and emission maxima at 528 and 617 nm, respectively.<sup>1</sup> Ethidium homodimer-1 has commonly been used to stain DNA in agarose gel electrophoresis and dead cells in cell viability assays.<sup>2-4</sup>

### References

1. Sabnis, R.W. *Handbook of biological dyes and stains: Synthesis and industrial applications*. 1st ed., John Wiley & Sons, Inc. Madison, NJ, USA (2010).
2. Zakharov, S.F., Garner, M.M., and Chrambach, A. Fluorescent labeling of DNA with ethidium homodimer without measurable decrease in DNA mobility: Application to automated gel electrophoresis apparatus. *Anal. Biochem.* **224**(1), 195-198 (1995).
3. Lévesque, A., Paquet, A., and Pagé, M. Measurement of tumor necrosis factor activity by flow cytometry. *Cytometry* **20**(2), 181-184 (1995).
4. Edwards, J.R., Diamantakos, E.A., Peuler, J.D., et al. A novel method for the evaluation of proximal tubule epithelial cellular necrosis in the intact rat kidney using ethidium homodimer. *BMC Physiol.* **7**, 1 (2007).

#### 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, 01/28/2026

#### 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