

# PRODUCT INFORMATION



## DAPI (hydrochloride) (solution)

Item No. 40796

**CAS Registry No.:** 28718-90-3  
**Formal Name:** 2-[4-(aminoiminomethyl)phenyl]-1H-indole-6-carboximidamide, dihydrochloride  
**Synonyms:** 4',6'-Diamidino-2-phenylindole, FxCycle Violet

**MF:** C<sub>16</sub>H<sub>15</sub>N<sub>5</sub> • 2HCl

**FW:** 350.3

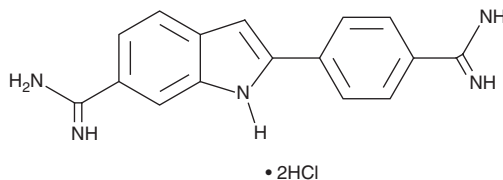
**Purity:** ≥98%

**Ex./Em. Max:** 347/453 nm

**Supplied as:** A solution in DMSO

**Storage:** -20°C

**Stability:** ≥2 years



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

### Description

DAPI is a fluorescent DNA probe.<sup>1,2</sup> It forms a fluorescent complex by attaching in the minor groove of adenine and thymine-rich sequences of DNA.<sup>1</sup> DAPI displays excitation/emission maxima of 347/453 nm, respectively. It has been commonly used as a cell viability dye and as a counterstain for DNA and chromosomes in fluorescent microscopy and flow cytometry applications.<sup>1,2</sup>

### References

1. Kapuscinski, J. DAPI: A DNA-specific fluorescent probe. *Biotech. Histochem.* **70(5)**, 220-233 (1995).
2. Zink, D., Sadoni, N., and Stelzer, E. Visualizing chromatin and chromosomes in living cells. *Methods* **29(1)**, 42-50 (2003).

#### 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, 03/19/2024

#### 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