

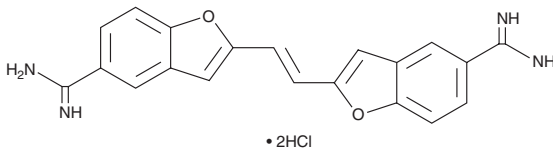
# PRODUCT INFORMATION



## True Blue

Item No. 26863

**CAS Registry No.:** 71431-30-6  
**Formal Name:** 2,2'-(1E)-1,2-ethenediylbis-5-benzofurancarboximidamide, dihydrochloride  
**Synonym:** NCI 240899  
**MF:** C<sub>20</sub>H<sub>16</sub>N<sub>4</sub>O<sub>2</sub> • 2HCl  
**FW:** 417.3  
**Purity:** ≥90%  
**Ex./Em. Max:** 373/404 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.

### Laboratory Procedures

True blue is supplied as a solid. A stock solution may be made by dissolving the true blue in the solvent of choice. True blue is soluble in the organic solvent methanol, which should be purged with an inert gas.

### Description

True blue is a fluorescent neuronal retrograde tracer that labels the nucleus, nucleolus, cell body, proximal dendrites, and axons of neurons.<sup>1</sup> It displays excitation/emission maxima of 373/404 nm, respectively.<sup>2</sup>

### References

1. Bentivoglio, M., Kuypers, H.G., Catsman-Berrevoets, C.E., *et al.* Fluorescent retrograde neuronal labeling in rat by means of substances binding specifically to adenine-thymine rich DNA. *Neurosci. Lett.* **12(2-3)**, 235-240 (1979).
2. Kasten, F.H. Table 2.3. *Fluorescent and Luminescent Probes for Biological Activity: A Practical Guide to Technology for Quantitative Real-Time Analysis* 33 (1999).

#### 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, 11/07/2022

#### 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](http://WWW.CAYMANCHEM.COM)