

PRODUCT INFORMATION



HBC 530

Item No. 40295

CAS Registry No.: 156840-13-0

Formal Name: 4-cyano- α -[[4-[(2-hydroxyethyl)methylamino]phenyl]methylene]-benzeneacetonitrile

MF: $C_{19}H_{17}N_3O$

FW: 303.4

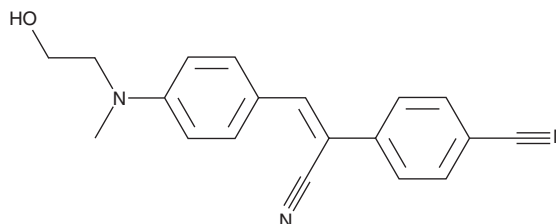
Purity: $\geq 98\%$

Ex./Em. Max: 485/530 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

HBC 530 is supplied as a solid. A stock solution may be made by dissolving the HBC 530 in the solvent of choice, which should be purged with an inert gas. HBC 530 is soluble (≥ 10 mg/ml) in DMSO.

Description

HBC 530 is a cognate dye ligand for the Pepper RNA aptamer.^{1,2} Upon binding to the Pepper RNA aptamer, the Pepper-HBC 530 complex displays excitation/emission maxima of 485/530 nm, respectively.¹ HBC 530 has been used in the live cell imaging of RNA.

References

1. Chen, X., Zhang, D., Su, N., *et al.* Visualizing RNA dynamics in live cells with bright and stable fluorescent RNAs. *Nat. Biotechnol.* **37**(11), 1287-1293 (2019).
2. Huang, K., Chen, X., Li, C., *et al.* Structure-based investigation of fluorogenic Pepper aptamer. *Nat. Chem. Biol.* **17**(12), 1289-1295 (2021).

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, 04/02/2024

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