

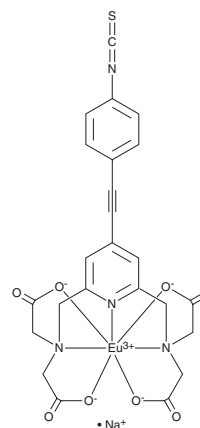
# PRODUCT INFORMATION



## Europium W1024 ITC

Item No. 31103

**CAS Registry No.:** 1204652-02-7  
**Formal Name:** [[N,N'-[[4-[[4-isothiocyanatophenyl]ethynyl]-2,6-pyridinediyl-κN]bis(methylene)]bis[N-[(carboxy-κO)methyl]glycinato-κN,κO]](4-)]-europate(1-), monosodium salt  
**Synonyms:** Eu-W1024 ITC, Europium W1024 Isothiocyanate  
**MF:** C<sub>24</sub>H<sub>18</sub>EuN<sub>4</sub>O<sub>8</sub>S • Na  
**FW:** 697.4  
**Ex./Em. Max:** 340/613 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Special Conditions:** Protect from light



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

### Description

Europium W1024 isothiocyanate (ITC) is a lanthanide-based fluorophore with a long emission lifetime and large Stokes shift.<sup>1,2</sup> It displays excitation/emission maxima of 340/613 nm, respectively.<sup>1</sup> Europium W1024 ITC contains a reactive isothiocyanate handle that reacts with primary amino groups at an alkaline pH. It is stable in the presence of high temperature, high EDTA concentration, or low pH but not in the presence of manganese. Europium W1024 ITC has commonly been conjugated to various molecules and used as a donor fluorophore in TR-FRET where its long emission lifetime provides a longer window of measurement for acceptor fluorescence and reduces the interference of autofluorescence.<sup>3-5</sup>

### Protein Labeling Protocol

The recommended conditions for labeling a protein are:

1. An amine-free buffer with a pH of 9.0-9.3
2. Overnight reaction at 4°C, protected from light
3. A molar excess of the europium chelate over protein to be labeled

*NOTE: The labeling efficiency will also depend on the protein concentration and its isoelectric point (pI). It is recommended that users perform a titration to identify the ideal chelate:protein molar ratios for their assay. A purification step is necessary to separate unreacted chelate from the labeled protein.*

It is recommended to store the chelate in an acidic buffer (pH 4-4.5) at a high concentration between -20 and -80°C.

### References

1. Mathis, G. and Bazin, H. *Stable luminescent chelates and macrocyclic compound*. Lanthanide luminescence: Photophysical, analytical and biological aspects. Hänninen, P. and Härmä, H, 1st, Springer, Berlin, Heidelberg (2010).
2. Mathis, G. *Clin. Chem.* **41(9)**, 1391-1397 (1995).
3. Moshinsky, D.J., Ruslim, L., Blake, R.A., et al. *J. Biomol. Screen.* **8(4)**, 447-452 (2003).
4. Scheepstra, M., Leysen, S., van Almen, G.C., et al. *Nat. Commun.* **6**, 8833 (2015).
5. Poore, D.D., Hofmann, G., Wolfe, L.A., 3rd., et al. *SLAS Discov.* **24(2)**, 175-189 (2019).

#### 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, 10/10/2025

#### 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