

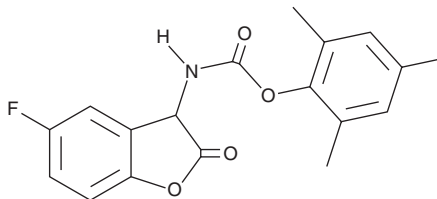
# PRODUCT INFORMATION



## 2-Coumaranone-1-L

Item No. 27389

**CAS Registry No.:** 1887057-34-2  
**Formal Name:** N-(5-fluoro-2,3-dihydro-2-oxo-3-benzofuranyl)-carbamic acid-2,4,6-trimethylphenyl ester  
**MF:** C<sub>18</sub>H<sub>16</sub>FNO<sub>4</sub>  
**FW:** 329.3  
**Purity:** ≥90%  
**Supplied as:** A crystalline 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

2-Coumaranone-1-L is supplied as a crystalline solid. A stock solution may be made by dissolving the 2-coumaranone-1-L in the solvent of choice, which should be purged with an inert gas. 2-Coumaranone-1-L is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of 2-coumaranone-1-L in these solvents is approximately 30 mg/ml.

2-Coumaranone-1-L is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2-coumaranone-1-L should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 2-Coumaranone-1-L has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

2-Coumaranone-1-L is a chemiluminescent probe.<sup>1</sup> In the presence of a base and oxygen, 2-coumaranone-1-L displays an emission maximum of 442 nm.

### Reference

- Schramm, S., Ciscato, L.F., Oesau, P., *et al.* Investigations on the synthesis and chemiluminescence of novel 2-coumaranones - II. *ARKIVOC* **2015(5)**, 44-59 (2015).

#### 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, 02/16/2024

#### 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