

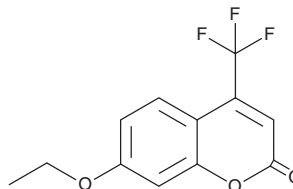
# PRODUCT INFORMATION



## 7-Ethoxy-4-(trifluoromethyl)coumarin

Item No. 34610

**CAS Registry No.:** 115453-82-2  
**Formal Name:** 7-ethoxy-4-(trifluoromethyl)-2H-1-benzopyran-2-one  
**Synonym:** EFC  
**MF:** C<sub>12</sub>H<sub>9</sub>F<sub>3</sub>O<sub>3</sub>  
**FW:** 258.2  
**Purity:** ≥95%  
**Ex./Em. Max:** 410/510 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

7-Ethoxy-4-(trifluoromethyl)coumarin is supplied as a solid. A stock solution may be made by dissolving the 7-ethoxy-4-(trifluoromethyl)coumarin in the solvent of choice, which should be purged with an inert gas. 7-Ethoxy-4-(trifluoromethyl)coumarin is slightly soluble in methanol and chloroform.

### Description

7-Ethoxy-5-(trifluoromethyl)coumarin is a fluorogenic substrate for cytochrome P450s (CYPs).<sup>1</sup> Upon O-deethylation by CYPs, 7-hydroxy-4-(trifluoromethyl)coumarin (HFC) is released and its fluorescence can be used to quantify CYP activity. HFC displays excitation/emission maxima of 410/510 nm, respectively.

### Reference

1. Buters, J.T.M., Schiller, C.D., and Chou, R.C. A highly sensitive tool for the assay of cytochrome P450 enzyme activity in rat, dog and man. Direct fluorescence monitoring of the deethylation of 7-ethoxy-4-trifluoromethylcoumarin. *Biochem. Pharmacol.* **46(9)**, 1577-1584 (1993).

#### 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, 10/17/2022

#### 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)