

## 7-amino-4-trifluoromethyl-coumarin

## Instruction Manual

Catalog Number	PK-CA577-1077-100
Description	The free form of AFC can be used as a fluorescent marker and standard in fluorometric caspase assays. Solublize the AFC powder in methanol or DMSO (1 mM) and then dilute it into aqueous buffer at different concentrations within the 0-4 $\mu$ M range. Read 100 $\mu$ l/sample.
Quantity	100 mg
Sequence / Molecular Weight / Molecular Formula	229,2 Da; C <sub>10</sub> H <sub>6</sub> NO <sub>2</sub> F <sub>3</sub>
Purity	>99% by TLC
Appearance / Formulation / Solubility	Yellow powder; soluble in methanol or DMSO (25 mg/ml).
Storage & Stability	Store at room temperature as supplied. After reconstitution, store at -20°C. Stable for one year.
Applications	NA
References	1. Fletcher, A., et al. (1977) Appl. Phys. 45:295-299. Bissel, E.R., et al. (1980) J. Org. Chem. 45:2283-2289.
Caution	NA

---

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.

---