

## Data Sheet

### *Fluorescein-labeled S-adenosyl homocysteine*

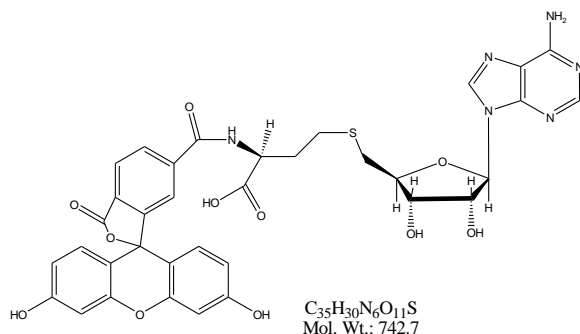
**Catalog : # 52012**

**Lot# : 160811**

**DESCRIPTION:** S-adenosyl homocysteine, fluorescein conjugate.

**PURITY:** >90% by HPLC

**STRUCTURE:**



**APPLICATIONS:** Study enzyme kinetics, and screen small molecular inhibitors of S-adenosyl homocysteine hydrolase and histone methyltransferases for drug discovery and HTS applications

**SUPPLIED AS:** DMSO solution

**QUANTITY:** 20  $\mu$ l @ 12.5  $\mu$ M

**STORAGE:** -20°C or -70°C

**WARNINGS:** Protect from light. Avoid freeze/thaw cycles.

**REFERENCES:**

1. Coward JK, Slisz EP. J Med Chem. 1973 May;16(5):460-3.
2. Finkelstein JD. Clin Chem Lab Med. 2007;45(12):1694-9.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)