



Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

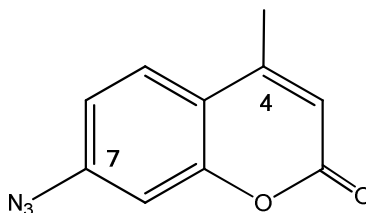
Technical Data Sheet

For research use only

Not intended or approved for
diagnostic or therapeutic use.

Product Name: 7-azido-4-methylcoumarin (AzMC)

Product Number: D-0010



Molecular Formula: C₁₀H₇N₃O₂ **MW:** 201.18

Solubility: Water, TRIS buffer

Storage: -20 °C

Field of Interest: This novel and selective probe for hydrogen sulfide (H₂S), 7-azido-4-methylcoumarin (AzMC), is useful for the monitoring of cystathionine β-synthase (CBS) catalyzed reactions, which plays a critical role in human sulfur metabolism. Deactivated CBS is characterized by high plasma levels of homocysteine and H₂S, resulting in clinical manifestations of numerous disease states. AzMC, used as a fluorogenic probe and concurrent CBS activity, can be used to select compounds for activation or inhibition of this enzyme, and is suitable for high-throughput screening and the development of potential therapeutics.

References: M.K. Thorson, T. Majtan, J.P. Kraus, A.M. Barrios "Identification of Cystathionine β-Synthase Inhibitors Using a Hydrogen Sulfide Selective Probe." *Angewandte Chemie (Int. ed. Eng.)*
DOI: 10.1002/anie.201300841

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. 7-Azido-4-methylcoumarin is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposures or risks.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.