

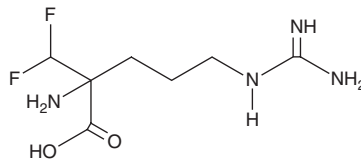
# PRODUCT INFORMATION



## $\alpha$ -(difluoromethyl)-DL-Arginine

Item No. 16415

**CAS Registry No.:** 69955-43-7  
**Formal Name:** 2-(difluoromethyl)-arginine  
**Synonyms:** DFMA, RMI 71897  
**MF:** C<sub>7</sub>H<sub>14</sub>F<sub>2</sub>N<sub>4</sub>O<sub>2</sub>  
**FW:** 224.2  
**Purity:**  $\geq 95\%$   
**UV/Vis.:**  $\lambda_{\max}$ : 202 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

$\alpha$ -(difluoromethyl)-DL-Arginine (DFMA) is supplied as a crystalline solid. Aqueous solutions of DFMA can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of DFMA in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Bacteria synthesize the cellular growth factor putrescine through a number of pathways. One pathway involves the decarboxylation of arginine by arginine decarboxylase to produce agmatine, which is then degraded to putrescine. While important for various cellular processes (e.g., cell division, differentiation, environmental stress responses) in all living organisms, including plants, high levels of this polyamine can be toxic.<sup>1</sup> DFMA is an enzyme-activated, irreversible inhibitor of the arginine decarboxylases of *E. coli* ( $K_i = 800 \mu\text{M}$ ), *P. aeruginosa*, and *K. pneumoniae*.<sup>2</sup> At 0.01 mM, it has been shown to prevent the osmotic stress-induced increase in arginine decarboxylase activity and putrescine synthesis in oat leaf cells.<sup>3,4</sup> DFMA can also reduce the growth of *T. cruzi* in mammalian host cells at a minimal concentration of 10 mM and prevent the growth of *C. parvum* in a T cell receptor alpha-deficient mouse model when combined with various polyamine analogs.<sup>5,6</sup>

### References

1. Gill, S.S. and Tuteja, N. Polyamines and abiotic stress tolerance in plants. *Plant Signal. Behav.* **5**(1), 26-33 (2010).
2. Kallio, A., McCann, P.P., and Bey, P. DL- $\alpha$ -(Difluoromethyl)arginine: A potent enzyme-activated irreversible inhibitor of bacterial decarboxylases. *Biochemistry* **20**(11), 3163-3168 (1981).
3. Tiburcio, A.F., Kaur-Sawhney, R., and Galston, A.W. Polyamine metabolism and osmotic stress. II. Improvement of oat protoplasts by an inhibitor of arginine decarboxylase. *Plant Physiol.* **82**(2), 375-378 (1986).
4. Flores, H.E. and Galston, A.W. Polyamines and plant stress: Activation of putrescine biosynthesis by osmotic shock. *Science* **217**(4566), 1259-1261 (1982).
5. Kierszenbaum, F., Wirth, J.J., McCann, P.P., et al. Arginine decarboxylase inhibitors reduce the capacity of *Trypanosoma cruzi* to infect and multiply in mammalian host cells. *Proc. Natl. Acad. Sci. USA* **84**(12), 4278-4282 (1987).
6. Yarlett, N., Waters, W.R., Harp, J.A., et al. Activities of DL- $\alpha$ -difluoromethylarginine and polyamine analogues against *Cryptosporidium parvum* infection in a T-cell receptor alpha-deficient mouse model. *Antimicrob. Agents Chemother.* **51**(4), 1234-1239 (2007).

#### 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, 11/14/2022

#### 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