



014PR-01

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Adenine Deaminase

A Recombinant Enzyme

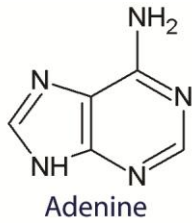
(Cat. # 786-428)



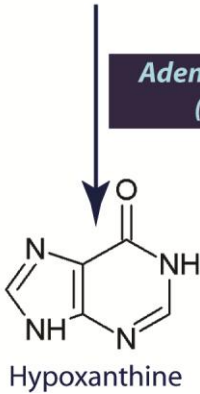
think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

Adenine Deaminase (EC 3.5.4.2) converts Adenine to Hypoxanthine. A highly purified recombinant enzyme protein supplied at a 2 μ M concentration.



Adenine



Hypoxanthine

Adenine Deaminase
(EC 3.5.4.2)

ITEM(S) SUPPLIED (Cat. # 786-428)

Description	Size
Adenine Deaminase (EC 3.5.4.2) [2 μ M]	0.25ml

STORAGE CONDITION

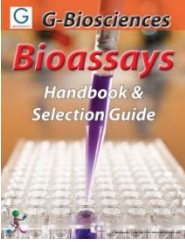
Shipped on dry ice. Upon receiving, store it at -80°C.

PROTOCOL

Reconstitute with pure water or a buffer of choice to desired concentration.

RELATED PRODUCTS

Download our Bioassays Handbook:



<http://info.gbiosciences.com/complete-bioassay-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 7/3/2014 CMH



www.GBiosciences.com