



570PR-01

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

A Geno Technology, Inc. (USA) brand name

Non Animal Protein Standards

(Cat. # 786-438, 786-439)



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

INTRODUCTION

G-Biosciences Non Animal Protein Standards are protein standards for use in protein estimations that are generated from non-animal proteins.

The prediluted Non Animal Protein Standards are ready to use solutions prepared in an aqueous buffer with a preservative for product stability. They are provided in six different concentrations ranging from 0.1mg/ml to 1.0mg/ml for running standard calibration in protein estimations.

ITEM(S) SUPPLIED

Cat. #	Description	Size
786-438	Non Animal Protein Standards [2mg/ml]	2 x 5ml
786-439	Non Animal Protein Standards (Prediluted) <i>Non Animal Protein Standards [0.1mg/ml]</i> <i>Non Animal Protein Standards [0.2mg/ml]</i> <i>Non Animal Protein Standards [0.3mg/ml]</i> <i>Non Animal Protein Standards [0.5mg/ml]</i> <i>Non Animal Protein Standards [0.8mg/ml]</i> <i>Non Animal Protein Standards [1.0mg/ml]</i>	5ml of each

STORAGE CONDITIONS

The kit is shipped at ambient temperature. Store refrigerated at 4°C, upon arrival. For long term storage keep at -20°C and is stable for 1 year. Avoid frequent freeze-thaw.

PROTOCOL

Non Animal Protein Standards (Prediluted)

For the prediluted Non Animal Protein Standards, use the appropriate volume direct in the protein assay. 50µl for 1ml protein assays and 5µl for microwell assays is recommended.

Non Animal Protein Standards [2mg/ml]

For minimizing interference, it is important to prepare the appropriate diluted protein standard in the same diluent used for the test protein sample.

For the Dilute Protocol, prepare a 0.1mg BSA/ml stock solution by mixing 50µl 2mg/ml stock with 950µl diluent. Use this stock for preparing diluted protein standard for the micro protocol assay.

For Standard Protocol (25-2000 μ g/ml)

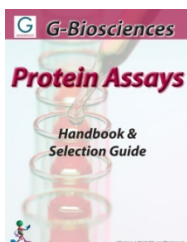
2mg/ml Standard (μ l)	Diluent (μ l)	Final Concentration (μ g/ml)
400	0	2000
300	100	1500
200	200	1000
150	250	750
100	300	500
50	350	250
25	375	125
5	395	25
0	400	0 (Blank)

For Dilute Protocol (2.5-25 μ g/ml)

0.1mg/ml Standard (μ l)	Diluent (μ l)	Final Concentration (μ g/ml)
250	750	25
200	800	20
150	850	15
100	900	10
50	950	5
25	975	2.5
0	1000	0 (Blank)

RELATED PRODUCTS

Download our Protein Assay Handbook



<http://info.gbiosciences.com/complete-protein-assay-guide/>

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



www.GBiosciences.com