



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ technical@GBiosciences.com

CERTIFICATE OF ANALYSIS

Product: RAPIDstain™

Size: 1L

Catalog Number: 786-31

Lot Number: 201812

TEST - Q.C. GEL STAINING

SDS-PAGE

Gradient 15 well Gel 4-20%, SDS-Tris-Glycine

Protein BSA

BSA Loaded/Well:

1 μ g, 0.5 μ g, 0.25 μ g, 0.125 μ g, 0.062 μ g, 0.031 μ g, 0.015 μ g (15ng), 0.0078 μ g (8ng), 0.0039 μ g (4ng)

BAND VISIBILITY TEST:

1 μ g - 0.031 μ g bands visible within 2-3 minutes	OK	Yes / No
1 μ g - 0.015 μ g bands visible within 5 minutes	OK	Yes / No
8ng - 4ng bands visible within 30 minutes	OK	Yes / No
Bands visibility after washing (1 μ g - 8ng)	8ng to 1 μ g protein bands visible	

This document has been produced electronically and is valid without signature

Date of Print: December 29, 2020



think proteins! think G-Biosciences!

www.GBiosciences.com