



Certificate of Analysis

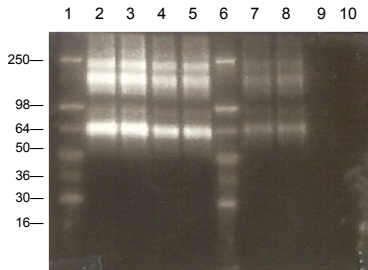
Product: Anti-Phosphotyrosine-HRP Mouse MAb V 1.0
Catalog #: APY03-HRP
Product Description: Mouse monoclonal IgG2b
Clone 27B10.4
Purified by protein G affinity chromatography
Amount: 100 µl per tube when reconstituted (APY03-HRP)
25 µl per tube when reconstituted (APY03-S-HRP)
Lot: 011

TEST	SPECIFICATON	LOT RESULTS
Appearance	White lyophilized powder	White lyophilized powder
Protein Quantitation Protein quantitated with Precision Red protein assay reagent	60-120 µg per tube Cat # APY03-HRP	80 µg
	15-30 µg per tube Cat# APY03-S-HRP	20 µg
Aliquot size is determined by biological activity therefore µg amounts per tube will vary between lots.		
Sensitivity/Specificity	10 ng of chemically labeled phosphotyrosine-BSA detected on western blot by chemiluminescence when 1:6000 Ab dilution and 30 second exposure time is used. A 1,000 ng sample of unlabeled BSA is not detected under identical conditions.	10 ng phosphotyrosine-BSA detected. 1,000 ng BSA not detected
Endogenous specificity	Detects multiple phosphoproteins in western blot analysis of 10 µg of cell lysate from NIH3T3 cells treated with pervanadate. Extremely faint or no signal is detected in untreated NIH3T3 or calyculin A treated A431 cell lysates. APY03 used at 1:6000 dilution, 10 second exposure for chemiluminescence detection.	Multiple phosphoproteins detected in pervanadate treated NIH3T3 cell lysates. Faint or no signal in untreated or calyculin A treated cell lysates.
QC Release Date	5/28/15	

Kim Middleton
QC Manager

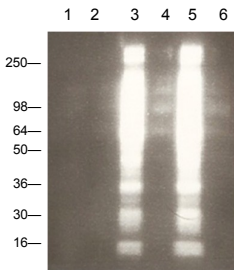
Lot specific QC Data

Sensitivity/Specificity: Lot 011



Lane 1 & 6: Markers, Lanes 2 -3: 10 ng phosphotyrosine labeled BSA, Lanes 4-5: 5 ng phosphotyrosine labeled BSA, Lanes 7-8: 1 ng phosphotyrosine labeled BSA, Lanes 9-10: 1000 ng unlabeled BSA. See data sheet for western blot protocol. A 1:6,000 dilution of APY03-HRP was used to develop the western blot. Bands were viewed by chemiluminescence, 30 second exposure time.

Endogenous Sensitivity/Specificity: Lot 011



Lane 1: 10 µg cell lysate from calyculin A treated A431 cells, Lane 2: 10 µg cell lysate from untreated A431 cells, Lanes 3 & 5: 10 µg and 20 µg respectively of cell lysate from pervanadate treated NIH3T3 cells (see data sheet for method), Lanes 4 & 6: 10 µg and 20 µg respectively of cell lysate from untreated NIH3T3 cells (see data sheet for method). A 1:6,000 dilution of APY03-HRP was used to develop the western blot. Bands were viewed by chemiluminescence, 10 second exposure time.