











PRODUCT DATA SHEET

Ready-to-Screen Tissue BLOTS™ ***Single Tissue Blot - Single Species***

Catalog #: TB68

Lot #: -----

Components: Protein blot of Normal Human Thyroid Single Tissue sample arranged as follows:

Lane 1	Protein Marker*	211,806		Myosin
Lane 2	Human Thyroid (Normal)	121,020		β-galactosidase
		100,216		Bovine Serum Albumin
		54,395		Ovalbumin
		38,708		Carbonic Anhydrase
		29,806		Soybean Trypsin Inhibitor
		20,040		Lysozyme
		7,331		Aprotinin

Size: 1 Blot

Storage Condition: 4° C

* Lot #: 300002325-BR

Methods Involved: The proteins were isolated from normal human thyroid tissue by preparing a tissue homogenate in the presence of protease inhibitors. Protein sample (50µg) from normal human thyroid tissue, solubilized in SDS-lysis buffer was electrophoresed in a 10 well, 4-20% SDS-polyacrylamide gradient gel, followed by electroblotting on PVDF membrane.

Quality Control: Proteins isolated from each lot were run on 4-20% gel and stained with G-Biosciences *RapidStain™* to check for its quality. Actin antibody was used to test the separation and transfer of protein from each lot.

Instructions for Use: Remove the blot from the pouch and wash with an appropriate buffer (1X TBST or PBST) 1-2 times. Block the membrane with a protein blocking agent; e.g., G-Biosciences *NAP™*-Blocker or *BLOT-QuickBlocker™*, and incubate with the primary and secondary antibodies diluted in blocking solution, following the standard protocol. Develop the blot with chemiluminescent or chromogenic detection reagents for the detection of the specific protein.

Rev 12.08.08-SA/MM/IA



think proteins! think G-Biosciences!

www.GBiosciences.com