

**D1000.IC Lot No. 2310034**

Beagle Dog Intestine Cytosol

Untreated, Male, Pool of 6

1.0 mL at 4 mg protein / mL

Suspension medium: 50 mM Tris·HCl, 150 mM KCl, 1 mM EDTA, 20% glycerol, heparin, PMSF, leupeptin, DTT, aprotinin

**Animal Information**

Species: Dog; *Canis familiaris*  
Strain: Beagle  
Sex: Male  
Age: >6 months  
Vendor: Marshall BioResources, North Rose, NY

Animals were housed in an AAALAC-accredited facility and allowed to acclimate  $\geq$  seven days before use.

Food: Nutrena (*ad libitum*)  
Water: Automatic watering system, tap water (*ad libitum*)  
Light/dark cycle: Not monitored  
Temperature: Ranges from 62°-82°F  
Humidity: Not monitored  
Cage: Indoor/outdoor run cages, plastic coated rod bottom, sanitized at least every 2 weeks

Aprotinin, Leupeptin, and Phenylmethylsulfonyl-fluoride were used in the preparation of this cytosol. Subcellular fractions were prepared from duodenal and jejunal tissue.

**Store at -80°C**

CAUTION: This sample should be considered as a potential biohazard and universal precautions should be followed. Intended for *in vitro* use only.

These data were generated by and are the property of Xenotech. These data are not to be reproduced, published or distributed without the express written consent of Xenotech.

This data sheet serves as a Certificate of Analysis and has been approved by **Stephanie Helmstetter, Assistant Director.**

Signature and Date: Stephanie Helmstetter 03 February 2023