

D1000.IC Lot No. 2110216

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: Envigo, Cumberland, VA

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.

Datasheet prepared 09 July 2021