# **DATA SHEET**





## CMV pp52 (UL44) 51 kDa

Cytomegalovirus pp52 (UL44) protein recombinant, *E. coli* 

Cat. No.	Amount
PR-1443-S	200 μg
PR-1443-L	1 mg

#### For general laboratory use.

**Shipping:** shipped on gel packs **Storage Conditions:** store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 6 months

Molecular Weight: about 51 kDa

Form: liquid (Supplied in Tris-HCl 50 mM pH 8.0, Sodium Chloride 60

mM, Glutathione 10 mM, Sarcosyl 0.25 %, Glycerol 50 %)

## **Applications:**

Suitable for use in ELISA. Each laboratory should determine an optimum working titer for use in its particular application. Other applications have not been tested but use in such assays should not necessarily be excluded.

#### **Description:**

The protein contains immunodominant region of CMV pp52 (UL44), amino acids 202-434. The protein is fused to a GST tag.

### Specificity:

Immunoreactive with sera of CMV-infected individuals