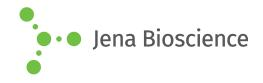
DATA SHEET





Varicella Zoster Virus (VZV) Glycoprotein Antigen

Strain Ellen native, HEL-299 cell culture

Cat. No.	Amount
PR-BA104VSG-S	250 μg
PR-BA104VSG-L	1 mg

For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at -80 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Form: liquid

Culture system: HEL-299 cell culture

Protein concentration: Lot specific. Determined by Bradford assay (Biorad) using BSA as standard.

assay (Biorau) using BSA as standard.

Preservation: None

Analysis: ELISA

Preparation: Infected cells are extracted by detergent treatment and the glycoproteins are isolated by Lectin affinity chromatography. The resulting antigen is presented in glycine buffer (pH 9.5). Residual amounts of detergent may be present.

Inactivation: This product has been treated with detergent. Confirmed inactive by attempted growth under original culture conditions. Since no procedure can guarantee absolute sterility, the reagent should be handled with appropriate safety precautions.