





Yersinia enterocolitica O:9 Antigen

native, liquid culture

Cat. No.	Amount	Culture system: Liquid culture
PR-BA13809-S	250 µg	Protein concentration: Lot specific. Determined by B assay (Biorad) using BSA as standard.
PR-BA13809-L	1 mg	
For general laboratory use.		Preservation: None
		Analysis: SDS-PAGE and ELISA
Shipping: shipped on dry ice		Preparation: Yersinia is grown in liquid culture and st to secrete YOPs (Yersinia outer proteins). YOP proteins cond from the culture medium are presented in phosphate buffere
Storage Conditions: store at -80 °C		
Additional Storag	e Conditions: avoid freeze/thaw cycles	
Shelf Life: 12 months		Inactivation: This product has been inactivated by s tration.Confirmed inactive by attempted growth under
Form: liquid (Supplied in PBS)		culture conditions. Since no procedure can guarantee sterility, the reagent should be handled with appropriat precautions.

