

## Product Data Sheet/Certificate of Analysis

### 0.5mL Zeba Column

**Compound:** 0.5mL Zeba Column  
**Storage:** Store at 4°C. Do Not Freeze



Catalog Number:	S-4024	Purity:	Nuclease Free; 50% gel slurry
Lot Number:		Expiration:	
Components:	High performance polymeric resin contained in a polypropylene column w/lid	Size:	

## Product Description

Zeba columns are used in the buffer exchange of proteins and desalting.

## Comments

Store at 4°C. Not for internal or external use in humans.

