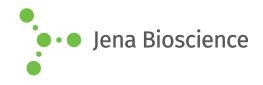
DATA SHEET





Agarose (blank)

| Cat. No. | Amount |
|----------|--------|
| AC-001S | 1 ml |
| AC-001L | 5 ml |

| | Agarose characteristics |
|------------------------------|---|
| Bead/Particle size | 45-165 μm |
| Recommended linear flow rate | 11.5 cm/h |
| Maximum pressure | 0.25 bar (3.6 psi) |
| pH stability | short term: 4 - 9 / long term: 4 - 9 |
| Chemical stability | Stable to all solutions commonly used in gel filtration including 8 M urea and 6 M guanidine hydrochloride Not stable in organic solvents! |
| Sterilization | Not autoclavable! |

Applications:

Suitable for control experiments with modified agarose.

Degree of substitution: -

Storage buffer: 20 % Ethanol

For general laboratory use.

Shipping: shipped at 4 °C

Storage Conditions: store at 4 $^{\circ}\text{C}$

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery