

EtOHNH₂-Agarose

Control gel

Ethanolamine-Agarose can be used as a control for unspecific adsorption effects.

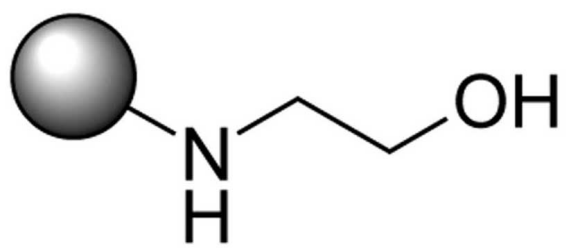
Ordering Information

[Order Online »](#)

| | |
|-------------|-------|
| BLG-E010-25 | 2.5ml |
| BLG-E010-60 | 6ml |

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|--------------------------|--|
| Use/Stability | Storage buffer should contain azide to prevent microbial growth. |
| Long Term Storage | +4°C |
| Shipping | Ambient Temperature |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|---------------------------------------|--|
| Alternative Name | Ethanolamine, immobilized on agarose gel |
| Couple Type | Modified nucleotides |
| Formulation | Storage buffer: 30mM phosphate, pH 7, in 20mg/l sodium azide. |
| Technical Info / Product Notes | For the Original Manufacturer's data sheet please click here . |



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com