

Heat Pretreatment Solution Citric

Ordering Information

[Order Online »](#)

| | |
|------------------|--------|
| ZTV-PT-0001-1000 | 1000ml |
|------------------|--------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|--------------------------|---|
| Use/Stability | Store, tightly closed, in the original vessel at temperatures between 2°C and 8°C |
| Long Term Storage | +4°C |
| Shipping | Blue Ice |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|---------------------------------------|--|
| Appearance | liquid; colorless |
| Formulation | Solution of salt in water |
| Technical Info / Product Notes | For the Manual please click here . |



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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