

# Diaphorase

Diaphorase belongs to a ubiquitous class of flavin-bound enzymes that catalyze the reduction of various dyes which act as hydrogen acceptors from the reduced form of di- and tri- phosphopyridine nucleotides (i.e. NADH, NADPH).

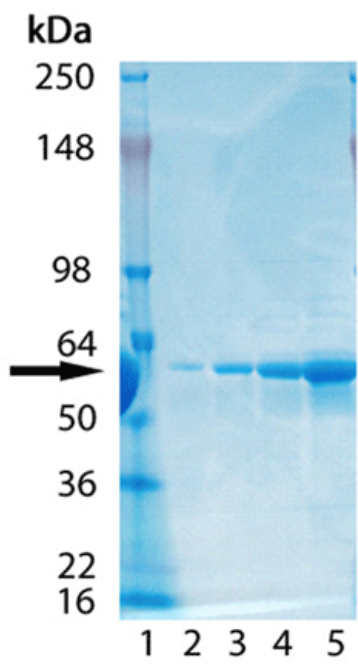
## Ordering Information

[Order Online »](#)

|                 |      |
|-----------------|------|
| ENZ-PRT155-0500 | 500U |
|-----------------|------|

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Short Term Storage -80°C

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

|                  |  |
|------------------|--|
| Alternative Name | NADPH dehydrogenase                                  |
| CAS              | 9001-68-7  |
| Formulation      | Liquid. In 50mM TRIS, pH 8.0, containing 150mM NaCl. |
| MW               | ~56 kDa  |
| Purity           | ≥95%   |
| Purity Detail    | Purified by multi-step chromatography.               |
| Source           | Produced in <i>E. coli</i> .                         |

Last modified: May 29, 2024



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)