

## Product Information

|                           |  |
|---------------------------|--|
| <b>Construct:</b>         | HMGB1 (2-215(end)-Avi-His)   |
| <b>Concentration:</b>     | 0.20 mg/ml   |
| <b>Species:</b>           | Human  |
| <b>Formulated In:</b>     | 8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol. |
| <b>Expression System:</b> | HEK293   |
| <b>Format:</b>            | Aqueous buffer solution  |
| <b>Stability:</b>         | At least 6 months at -80°C. Avoid freeze/thaw cycles.              |
| <b>Storage:</b>           | -80°C  |
| <b>Genbank Accession:</b> | NM_002128  |
| <b>MW:</b>                | 25 kDa   |
| <b>Glycosylation</b>      | This protein runs at higher M.W. by SDS-PAGE due to glycosylation. |
| <b>Purity:</b>            | ≥90%   |

## Quality Control Data

### 4-20% SDS-Page Coomassie Staining

