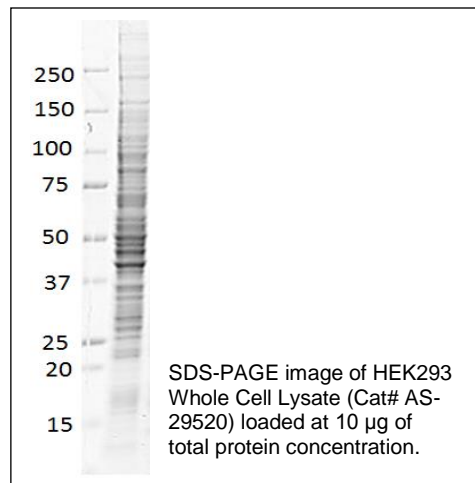


Update: September, 2017

## Product Data Sheet

---

|                    |   |
|--------------------|---|
| Product Name:      | HEK293 Whole Cell Lysate  |
| Catalog Number:    | AS-29520  |
| Lot Number:        | See label on vial   |
| Formulation:       | Lysate is in SDS-PAGE sample buffer (pH 6.8), heated for 5 minutes at 95-100 °C. The buffer is containing 62.5 mM Tris-HCl, 2% SDS, 10% glycerol, 0.002% bromophenol blue and reducing agent. |
| Size:              | 200 µg  |
| Concentration:     | 2 mg/mL   |
| Application Notes: | HEK293 Whole Cell Lysate is suitable for SDS-PAGE and/or Western blot. 10-30 µg of the total protein per lane is recommended.   |



**Storage:** Store at -80°C. Avoid repeating freeze-thaw cycles.

This product is for *in vitro* research use only.