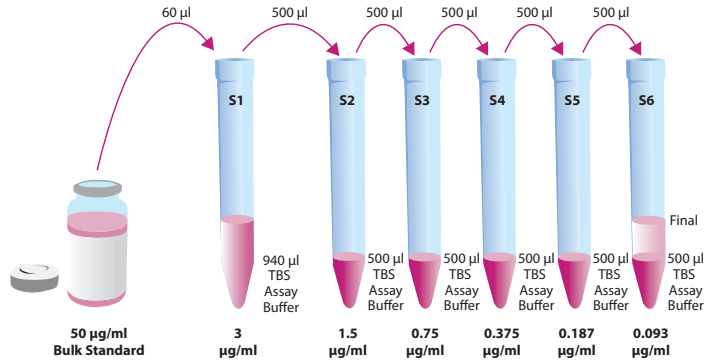


## His-Express Detection EIA Short Protocol

- TBS Buffer** - Dilute with 90 ml of UltraPure water.
- DEA Buffer** - Dilute the 2.5 ml vial to 25 ml or dilute the 12.5 ml vial to 125 ml of UltraPure water.
- AP Buffer** - Dilute the 5 ml vial to 750 ml or dilute the 12.5 ml vial to 1,875 ml of UltraPure water.
- Tracer** - Reconstitute the 100 dtn vial with 6 ml of TBS Buffer or the 500 dtn vial with 30 ml TBS Buffer.
- Antibody** - Reconstitute the 100 dtn vial with 6 ml of TBS Buffer or the 500 dtn vial with 30 ml TBS Buffer.
- Standard** - Prepare as described in the figure below.



| Steps                  | Reagent  | Blank  | NSB    | B <sub>0</sub> | Std/Sample |
|------------------------|--|--------|--------|----------------|------------|
| <b>1. Add Reagents</b> | TBS Buffer   | --     | 100 µl | 50 µl          | --         |
|                        | Standard/Sample  | --     | --     | --             | 50 µl      |
|                        | Tracer   | --     | 50 µl  | 50 µl          | 50 µl      |
|                        | Antibody   | --     | --     | 50 µl          | 50 µl      |
| <b>2. Incubate</b>     | <b>Cover plate and incubate 90 minutes at room temperature (22°C) on an orbital shaker</b> |        |        |                |            |
| <b>3. Wash</b>         | <b>Wash all wells four times</b>   |        |        |                |            |
| <b>4. Add Reagents</b> | pNPP Substrate Solution  | 200 µl | 200 µl | 200 µl         | 200 µl     |
| <b>5. Incubate</b>     | <b>Cover plate and incubate 30-60 minutes at room temperature</b>                          |        |        |                |            |
| <b>6. Read</b>         | <b>Read plate at a wavelength between 405-420 nm</b>                                       |        |        |                |            |

