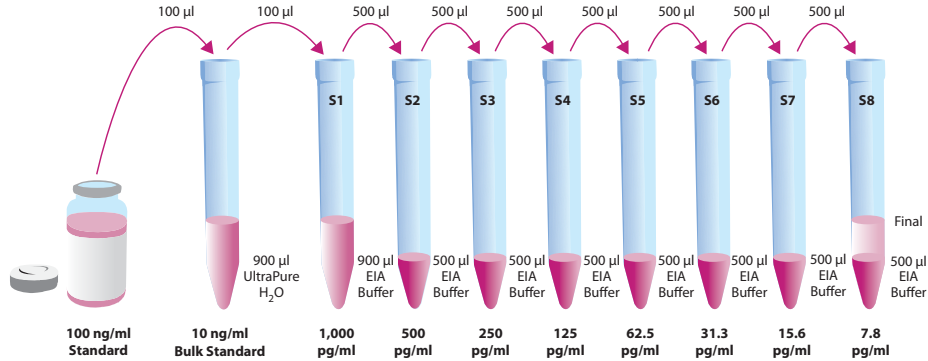


## Progesterone EIA Short Protocol

- EIA Buffer** - Dilute with 90 ml of UltraPure water.
- Wash Buffer** - Dilute the 5 ml vial to 2 L and add 1 ml Polysorbate 20 or dilute the 12.5 ml vial to 5 L and add 2.5 ml Polysorbate 20.
- Tracer** - Reconstitute the 100 dtn vial with 6 ml of EIA Buffer or the 500 dtn vial with 30 ml EIA Buffer.
- Antiserum** - Reconstitute the 100 dtn vial with 6 ml of EIA Buffer or the 500 dtn vial with 30 ml EIA Buffer.
- Standard** - Prepare as described in the figure below.



Steps	Reagent	Blank	TA	NSB	B <sub>0</sub>	Std/Sample
<b>1. Add Reagents</b>	EIA Buffer	--	--	100 µl	50 µl	--
	Standard/Sample	--	--	--	--	50 µl
	Tracer	--	--	50 µl	50 µl	50 µl
	Antiserum	--	--	--	50 µl	50 µl
<b>2. Incubate</b>	<b>Cover plate and incubate one hour at room temperature (22°C) with gentle shaking</b>					
<b>3. Wash</b>	<b>Wash all wells five times</b>					
<b>4. Add Reagents</b>	Tracer	--	5 µl	--	--	--
	Ellman's	200 µl	200 µl	200 µl	200 µl	200 µl
<b>5. Incubate</b>	<b>Cover plate and incubate 60-90 minutes RT with gentle shaking</b>					
<b>6. Read</b>	<b>Read plate at a wavelength between 405-420 nm</b>					

