

**PPAR $\alpha$ ,  $\delta$ ,  $\gamma$  Complete Transcription Factor Assay Short Protocol**

Steps	Reagent	Blk	NSB	PC	C1	S1-S44
1. Add reagents	CTFB	100 $\mu$ l	100 $\mu$ l	90 $\mu$ l	80 $\mu$ l	90 $\mu$ l
	Competitor dsDNA				10 $\mu$ l	
	Positive Control			10 $\mu$ l	10 $\mu$ l	
	Samples					10 $\mu$ l
2. Incubate	Cover plate and incubate overnight at 4°C without agitation					
3. Wash	Wash all wells five times					
4. Add reagents	Primary Antibody		100 $\mu$ l	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l
5. Incubate	Cover plate and incubate one hour at room temperature without agitation					
6. Wash	Wash all wells five times					
7. Add reagents	Goat Anti-Rabbit HRP conjugate		100 $\mu$ l	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l
8. Incubate	Cover plate and incubate one hour at room temperature without agitation					
9. Wash	Wash all wells five times					
10. Add reagents	Developer Solution	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l
11. Incubate	Monitor development in wells					
12. Add reagents	Stop Solution	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l
13. Read	Read plate at wavelength of 450 nm					

