



ARBOR
ASSAYS

Interactive Assay Solutions™

Prostaglandin E₂ (PGE₂) Multi-format Kits K051-H1/H5

As of June 1st, we no longer offer the K018-H, K018-HX or K018-C kits.

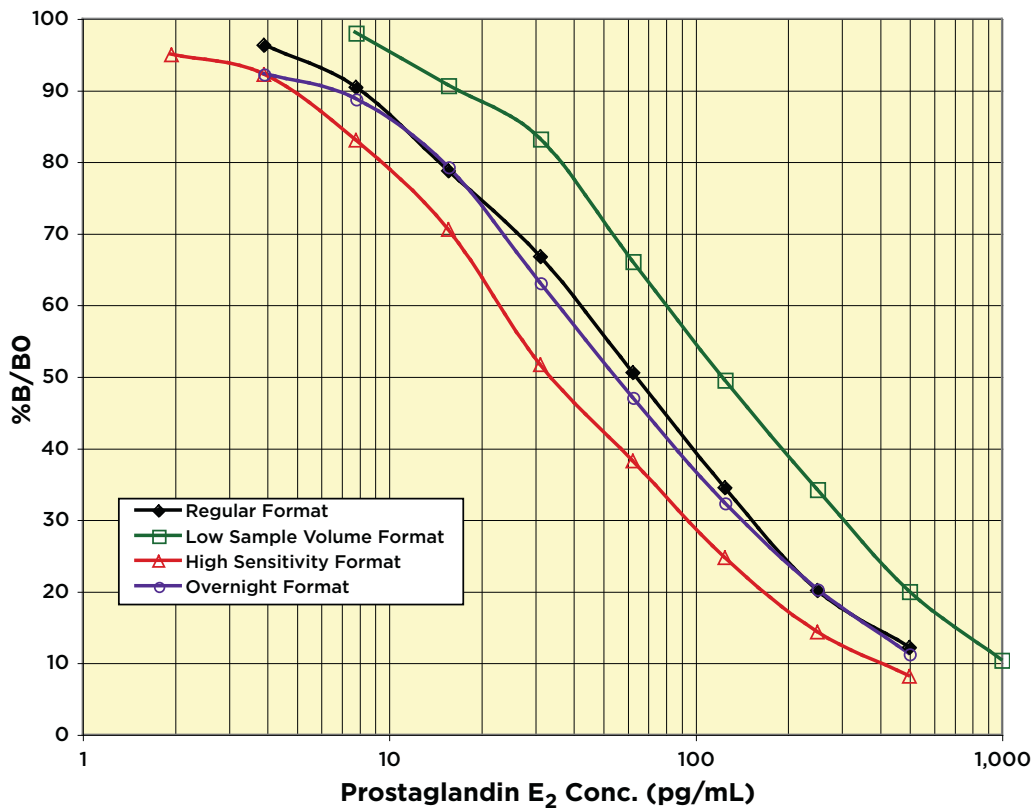
To help you become familiar with our new PGE₂ kit, we are extending our PGE₂ customers a free K051-H1 kit with your next order! Please let us know if you would like to try this kit for comparison purposes.

Our new PGE₂ EIA kit is radically different. The recently introduced DetectX® PGE₂ EIA kit, K051-H1 and K051-H5, uses an internally developed monoclonal to PGE₂. The antibody, which drives the performance of the assay has remarkable properties. The assay can handle all the formats and sensitivities of our K018-H, K018-HX and K018-C kits. Please see comparisons in the table below.

- Because of the rapid kinetics of the monoclonal antibody binding to PGE₂ the standard curve does not change if incubated for 2 hours at room temperature or overnight at 4°C.
- By increasing the volume of diluted sample into the well from 25 µL to 50 µL or 100 µL the sensitivity of the assay can be changed to cover all PGE₂ samples.
- The antibody has enhanced (lower) cross reactivity to PGE₁, PGF_{2α} and TXB₂.
- Mouse serum and plasma do not need to be extracted.

Samples are treated identically to the K018 kits. For the customer, this new kit gives them the knowledge that the one kit they purchase will be able to handle a very wide variety of sample needs. They can just increase the diluted sample volume from 25 µL to 50 µL or to 100 µL to get added sensitivity. If their lab procedure requires a 2 hour or an overnight assay the new K051-H kit will handle this change. It is also a colorimetric EIA with a reading at 450 nm, meaning that no special plate reader is needed.

We thank you for understanding, and feel that the K051-H kit will exceed your expectations!



Assay	Range (pg/mL)	Sensitivity (pg/mL)	Time to Answer	PGE ₁ Reactivity	PGF _{2a} Reactivity	TXB ₂ Reactivity
K051-H	1,000-1.95	3.1	2.5 hr or ON	27.3%	0.33%	<0.02%
K018-H	1,000-31.25	29.1	2.5 hr	108.9%	2.0%	0.3%
K018-HX	400-12.5	10.9	ON	108.9%	2.0%	0.3%
K018-C	320-5	4.8	ON	108.9%	2.0%	0.3%