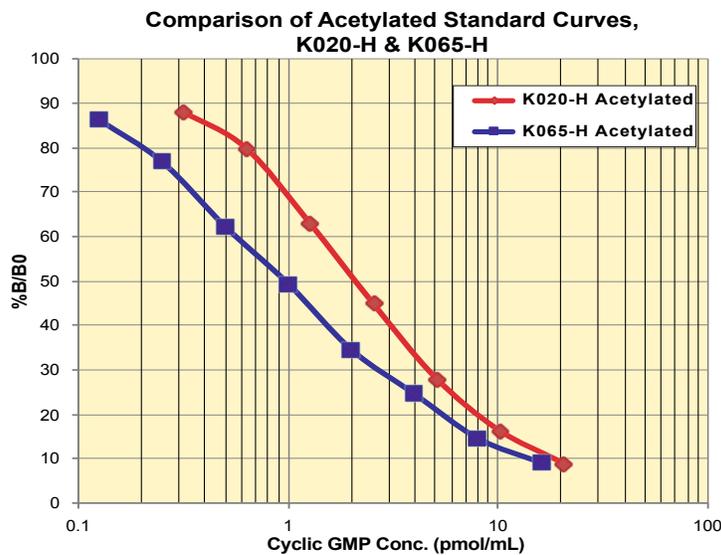




ABOUT THE KIT

When we invented Direct Cyclic Nucleotide EIA assays more than 20 years ago we knew they would be popular. The direct nature of the format for the assays allows cAMP & cGMP in cells to be measured in one step. At Arbor Assays we have refined the assay developing both colorimetric and chemiluminescent versions for cAMP and cGMP.

Cyclic GMP levels in most systems are an order of magnitude lower than those found for cyclic AMP and in our constant effort to deliver the easiest, most reliable and sensitive assays we have developed a new cGMP antibody to give you enhanced sensitivity. Our new Cyclic GMP EIA Kits (K065-H) are 2-fold more sensitive than our current cGMP assay (K020-H).



The improvement in sensitivity is shown by the Limits of Detection and calculated Sensitivity for the 2 assays, with K065-H being half that of K020-H.

Urine dilutions were run in the nonacetylated format of K065-H and K020-H, side by side, and measured cGMP concentrations were similar. Plasma dilutions were run in the acetylated format for each kit with only a slightly larger difference seen between the measured concentrations.

