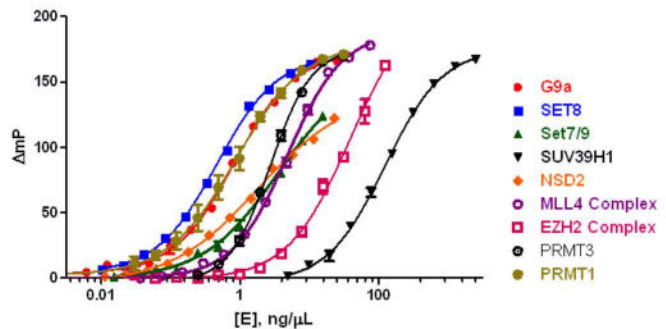


Ten Reaction Biology methyltransferases now validated with the Transcreener EPIGEN Methyltransferase Assay

BellBrook Labs and Reaction Biology are working together to make it easier to screen and profile methyltransferases. Reaction Biology has produced more than 30 recombinant human methyltransferases to the most stringent purity and activity requirements in the industry. We developed the Transcreener EPIGEN MT Assay to have the sensitivity needed to detect any MT at physiological SAM concentrations, and the robustness required for industrial HTS. And now, for each of Reaction Biology's enzymes, we are determining the optimal reaction conditions to produce an outstanding assay window under initial velocity conditions. This means you don't have to do any time consuming assay development. We will tell you how much enzyme to use, how long to run your reaction, and the right mix of detection reagents to get an assay window suitable for HTS. And the EPIGEN MT Assay has been validated on all of the major multimode readers, so we will tell you which filters and settings are right for your instrument. We have validated ten Reaction Biology enzymes thus far, and we are working hard on the rest so that you can spend less time on assay development and more on screening and profiling methyltransferases.

- [Learn more](#) about how BellBrook is validating Reaction Biology's methyltransferase enzymes to eliminate the need for assay development.
- [Visit](#) the Transcreener EPIGEN MT Assay webpage.
- See the recent [JBS paper](#) describing development and validation of the Transcreener EPIGEN Methyltransferase Assay.



For more information call toll free (866-313-7881) or email info@bellbrooklabs.com.