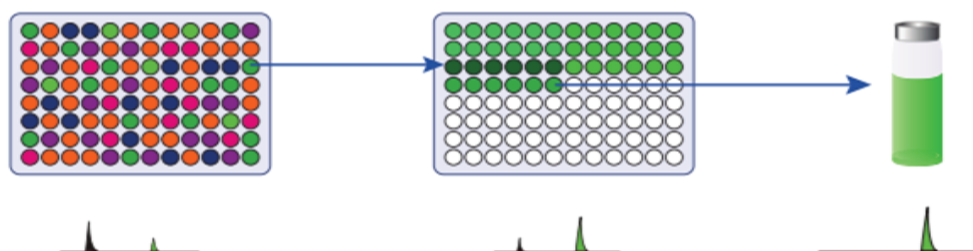




Metabolite Synthesis Made Easy

Identify & Produce Metabolites of Interest Quickly and Cost-Effectively



BMO Kit
Metabolite Production
Made Easy

Screen

Perform the primary screen. Select desired metabolite wells. Order the corresponding optimization kit.

Optimize

Perform the optimization kit. Identify the best production conditions. Order corresponding production kit.

Produce

Scale-up and isolate your metabolite

Metabolite

milligram quantities of metabolite generated

Product ID	Description	Product ID	Description	Product ID	Description
XTHCK1001-01	BMO Screening Kit	XTHCK1001-02	BMO Optimization Kit	XTHCK1001-03	BMO Production Kit

Features

- Utilizes state-of-the-art biomimetic technology, mimicking oxidative enzymes such as P450s, creating a synthetic chemical liver
- Innovative panels of catalytic chemical reaction conditions utilize organometallic catalysts in a multi-well, parallel format
- 3 simple steps to metabolite production
- Fast, cost-effective metabolite production in milligram quantities

Benefits

- Eliminate complex and costly synthetic chemistries
- Produce drug metabolites in a few days
- Rapid, inexpensive scale-up for NMR and *in vitro* testing
- Easy to implement - compatible with existing workflow

Contact us to learn more about our Metabolite Synthesis capabilities at info@xenotechllc.com or call us at **913.438.7450**.