



CryostaX® is XenoTech's single-freeze pooled human hepatocyte product, now available in plateable pools

*Gain more predictable and accurate results from pooled hepatocytes with CryostaX® products. XenoTech's unique approach measures every pool's characterization before and after pooling, for a double dose of quality control. Our patented single-freeze process provides the highest enzymatic activity possible by minimizing the additional cryoinjury caused by a second freeze required by conventional pooling methods.*

Pooled human hepatocytes are a preferred test system in many drug discovery and development applications. Like other pooled human hepatocytes, CryostaX® provides you with a representative population sample for predictable results; but our innovative process provides more extensive characterization, greater enzymatic activity and on-demand custom pooling with no minimum orders (does not include the plateable pools product). CryostaX® is available in 10- and 20- donor pools ideal for metabolism studies as well as 5-donor **plateable pools** ideal for long-term metabolism, drug transport and enzyme induction screening studies.

## Available CryostaX® Products:

### 10-Donor Pool: (HPCH10)

- 10 x 100µL = 1.0mL total volume
- AMY: 5 million cells

### 20-Donor Pool: (HPCH20-50)

- 20 x 50µL = 1.0mL total volume
- AMY: 5 million cells

### CryostaX® Plateable Pools (HPCH05+ & HPCH10+)

- 5-Donor Pool: 1.0mL volume; 5 million AMY
- 10-Donor Pool: 1.0mL volume; 5 million AMY

## What makes CryostaX® different?

### The single freeze difference

Hepatocytes are damaged by the freezing and thawing process. Sekisui XenoTech's unique, patented process used to prepare CryostaX®, eliminates the need for the second freeze used in conventional pooling methods, minimizing cryoinjury and preserving enzymatic activity. This assures you the highest quality available for dependable results.

## Free, on-demand, custom pooling, no minimum order

Design the CYP panel you need with precision. CryostaX® is flexible to your needs with no minimum order. XenoTech provides comprehensive, predictable outcomes for your custom pool, taking the guesswork out of your results. Design your own 10 or 20 donor custom pool for metabolism or induction purposes. (CryostaX® plateable pools)

## The most extensive characterization on the market

We stand by the quality of our products. We measure our enzymatic activities using validated LC/MS/MS methods and all characterization data are reviewed following a strict QC procedure. Each order of CryostaX® is accompanied by a lot-specific datasheet. We calculate theoretical enzymatic activities from the characterization of individual donors, and achieve those activities in the finished pooled product. All CYPs are targeted for average activity levels (based on our human hepatocyte characterization database) to assure lot-to-lot consistency.

## Characterization provided

CYP1A2	Phenacetin <i>O</i> -dealkylation
CYP2A6	Coumarin 7-hydroxylation
CYP2B6	Bupropion hydroxylation
CYP2C8	Amodiaquine <i>N</i> -dealkylation
CYP2C9	Diclofenac 4'-hydroxylation
CYP2C19	S-Mephenytoin 4'-hydroxylation
CYP2D6	Dextromethorphan <i>O</i> -demethylation
CYP2E1	Chlorzoxazone 6-hydroxylation
CYP3A4/5	Testosterone $\beta$ -hydroxylation
CYP3A4	Midazolam 1'-hydroxylation
UGT	7-Hydroxycoumarin glucuronidation
SULT	7-Hydroxycoumarin sulfonation



## CryostaX® Plateable Pools



Contact one of our Product Sales Specialists at [913.438.7450](tel:913.438.7450) or email them if you have questions or would like additional information on CryostaX®:

[info@xenotechllc.com](mailto:info@xenotechllc.com)

### Packaging

5-Donor Pool: (HPCH05+)

- 10 x 100µL = 1.0mL total volume
- AMY: 5 million cells

10-Donor Pool: (HPCH10+)

- 10 x 100µL = 1.0mL total volume
- AMY: 5 million cells

### Plateable Pools ideal for:

Long-term drug metabolism studies  
Drug transporter studies  
Enzyme Induction screening

### The CryostaX® Advantage

CryostaX® isn't a new procedure to add to your lab routine; it's a smarter and better way of optimizing your study results.

Preparation is as easy as pouring CryostaX® pellets into warmed media.

Your study deserves the best -- get better results with CryostaX®.  
Visit [www.xenotech.com](http://www.xenotech.com) or call **+913.438.7450** / in North America call toll-free **877.588.7530**