

XpressWay™ Profiles

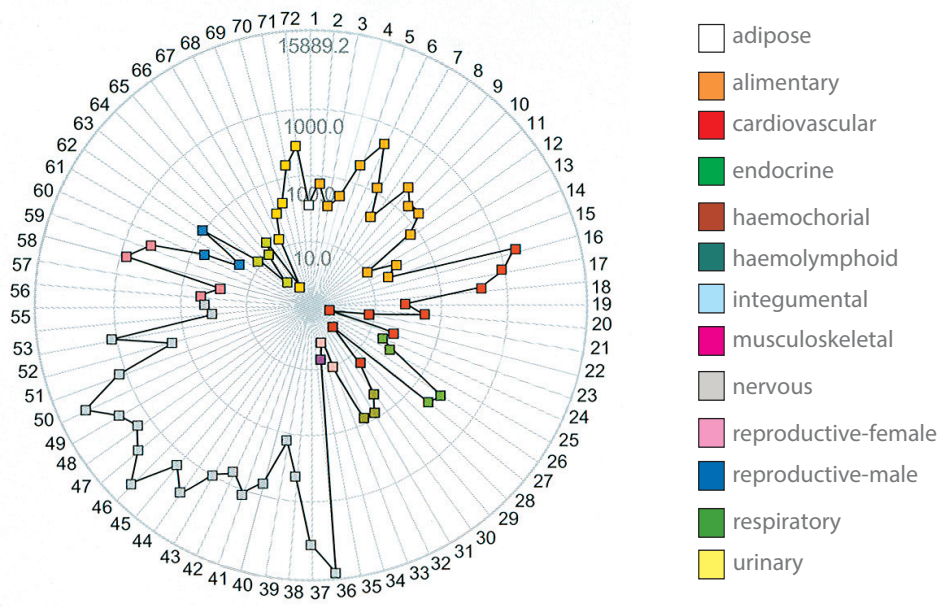
Snap-shot gene expression across major human organs and tissues

OVERVIEW

Asterand Bioscience's XpressWay™ Profiles offer quantitative gene expression data for a given target across 72 different non-diseased human tissue types, each from 3 different donors. The profiles provide a "snapshot" of where and to what extent a gene is expressed across the human body.

QUICK VISUAL IDENTIFICATION

XpressWay™ Profile data is presented as a stellar plot, enabling a fast and visual identification of where a given target is expressed. The 72 tissues are represented by the spokes of the wheel and the mean (n=3 donors/tissue) quantitative expression data are presented on a log scale from the center to the circumference of the wheel. The example below shows at a glance that GPR22 is most highly expressed in CNS and cardiovascular tissues.



HOW ARE XpressWay™ PROFILES GENERATED ?

- Xpressway™ Profiles are generated by qRT-PCR
- Asterand Bioscience designs all primer-probe sets and sequences are subject to homology searchers against the GenBank® database to confirm specificity for the target sequence
- Total RNA is extracted from each of the 72 tissue types (3 donors each), DNase treated and subject to the following QC:
 - Presence of 18S rRNA
 - Minimum copy numbers for control gene mRNA transcripts (qRT-PCR)
 - B-actin (amplicon length 295bp)- >3800 copies/100 ng RNA
 - GAPDH (amplicon length 751 bp) > 10,000 copies /100ng RNA
 - No DNA contamination

BENEFITS

- Excellent coverage of drugable targets and immediate access to data
- Comparison of human gene expression data across 72 normal tissue types
- Ability to access challenging targets and variants
- Comprehensive report includes donor demographics and clinical history, pathology, primer-probe sequences and all supporting interpretation of gene expression data
- Confirm therapeutic approach and exploit opportunities to switch indication

Issue Code: AB-PZ-XpressWay-1017

ABOUT ASTERAND BIOSCIENCE®

Asterand Bio is the leading global provider of high-quality, well-characterized human tissue and human tissue-based research solutions to drug discovery scientists. Our mission is to provide human-based solutions to accelerate the identification and validation of drug targets and enhance the selection of drug candidates with an increased likelihood of clinical success.

ETHICAL AND REGULATORY CONTROL

All human tissue samples are obtained with informed donor consent and are de-identified prior to banking and distribution. Asterand Bio acts in strict compliance with the US Health Insurance Portability and Accountability Act (HIPAA) and the United Kingdom Human Tissue Authority (HTA) and is accredited by the College of American Pathologists (CAP) and ISO 9001. Asterand Bio has GLP accreditation according to the requirements of Good Laboratory Practice (GLP)– Directive of the European Commission 2004/9/EEC, and a Good Clinical Laboratory Practice (GCLP) laboratory for clinical assay services.

RESEARCH SERVICES

- Immunohistochemistry
- ISH & FISH
- TMA, FDA & Custom
- Digital Pathology
- Histopathology
- qRT-PCR
- LCM
- siRNA Knockdown
- IHC/ISH Multiplexing
- Western Blotting
- Primary Cell Isolation
- 2D and 3D Culture
- Cell Based Assays
- Assay Development
- OrganDOT™
- Mutation Analysis

BIOMATERIALS AVAILABLE

- Fresh, Frozen & FFPE
- Normal & Diseased Tissue
- OCT Embedded
- Serum, Plasma & Blood
- Synovial Fluid
- RNA, DNA & Protein
- TMA, FFPE & Frozen
- Primary Cells
- Proprietary Cell Lines

THERAPEUTIC AREAS

- Oncology
- Cardiovascular
- Metabolic
- CNS
- Respiratory
- Inflammation
- Genitourinary
- Fibrosis

XpressBANK™ BIOREPOSITORY

The Asterand Bio XpressBANK™ Biorepository is an invaluable source of high-quality, well-characterized and validated tissue, biofluids, cell lines and molecular derivatives, all intended for research use in the drug and diagnostic development process. Our extensive repository contains over 200,000 human tissues and biofluids. This bank includes material that is prospectively obtained in a variety of formats including fresh frozen and FFPE, supported by comprehensive clinical data.

BioSPOKE™ CUSTOM BIOSPECIMEN PROCUREMENT

Through our BioSPOKE™ service, human tissues and clinical data are custom collected to meet unique research requirements within a wide range of disease categories or as normal controls. Samples are provided fresh or in multiple preserved formats. Specialized processing may be tailored to each research request. This service allows Asterand Bio to deliver human tissues, biofluids and accompanying clinical information to our clients that are collected in a manner which is most compatible with their experimental design.

PhaseZERO® RESEARCH SERVICES

For over 20 years, Asterand Bio's trusted scientists have worked collaboratively with clients to facilitate the efficient selection of drug targets and candidate therapeutics with the highest likelihood for clinical success by providing a variety of pre-clinical drug discovery services. From target and biomarker validation, characterization of therapeutic candidates for potency, effect and safety to supporting clinical diagnostic assay development and testing, Asterand Bio provides clients with high-quality experimental design, execution, data interpretation and reporting. Each project is managed by a dedicated scientific project manager and customized to meet each client's specific needs.



Unless otherwise noted, all trademarks are the property of Asterand Bioscience, Inc.
©2017 Asterand Bioscience, Inc. All rights reserved.

Please contact our Business Development Directors in the following regions to discuss your requirements:

United States

Tech One, Suite 501 | 440 Burroughs | Detroit, MI 48202 USA

T: 313-263-0960 | F: 313-263-0961

advantage@asterandbio.com

www.asterandbio.com

International

2A Orchard Rd. | Royston | Hertfordshire, SG8 5HD UK

T: +44 (0) 1763 211600 | F: +44 (0) 1763 211555



A BioIVT Company