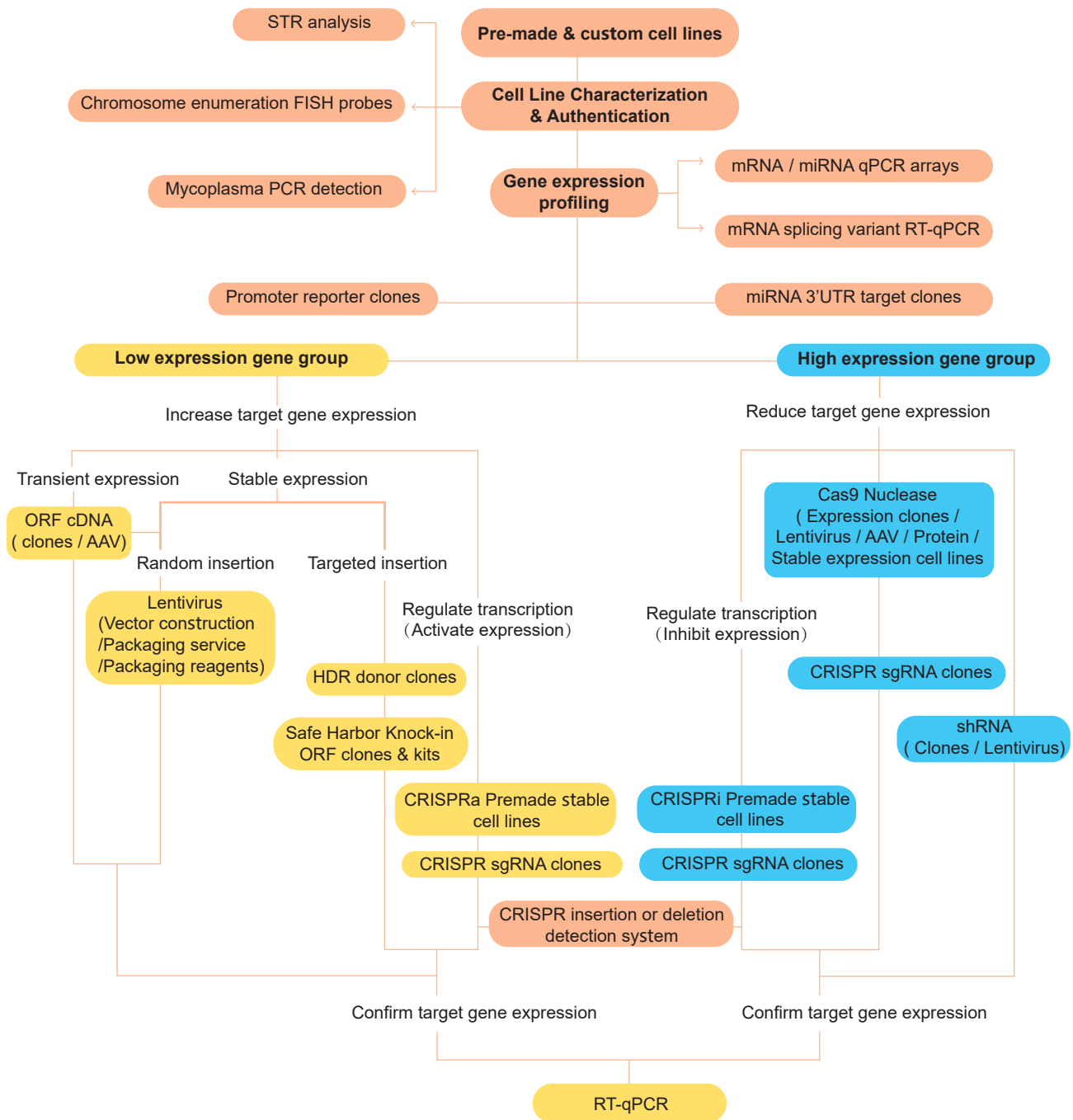


Functional Genomics - Companion Product Guide

○ Clickable link to products & services ○



♦ **Related Products & Services:** [Transfection Reagents] [AAV Particles] [Luciferase Assays] [Stable Cell Line Services]

♦ **More than 100,000 Publications**

Contact us:

GeneCopoeia, Inc.

9620 Medical Center Dr. Suite 101,
Rockville, MD 20850 USA

Email

inquiry@genecopoeia.com

Tel

+1 (301) 762-0888

Fax

+1 (301) 762-3888

Website

www.genecopoeia.com