

Product Information

Description:	URAT1 Cellular Transport Medium is an assay medium optimized for use with URAT1 HEK293 Cell Line in URAT1-based cellular transport assays.
Supplied As:	Aqueous solution.
Applications:	Use in URAT1 fluorescent transport assays with URAT1 HEK293 Cell Line.
Storage/Stability:	Upon receipt, store at 4°C. If unopened, it is stable for 6 months from date of receipt when stored as directed. Do not freeze!
Instructions for Use:	This is a ready-to-use buffer.
Optimized for use with:	URAT1 HEK293 Cell Line (BPS Bioscience #84223).