

ESTABLISHMENT AND CHARACTERIZATION OF A KIDNEY-DRUG INTERACTION MODEL BY STABLY EXPRESSING HOAT1 IN HEK 293T/17 CELLS

ATCC

Aaron Briley, BS, James Clinton, PhD, Brian Shapiro, PhD, Virginia Takahashi, MS, and Chaozhong Zou, PhD

ABSTRACT

Clearance of organic toxins by the kidney is a critical mechanism for mammalian homeostasis and for testing the toxicity of experimental drugs and other compounds. In this study, we created a HEK293T/17 cell line that stably expresses the human organic anion transporter (OAT1) and tested its transport capabilities.

INTRODUCTION

In vivo studies have shown that kidney membrane transporters play a key part in drug disposition and renal clearance. One such transporter is OAT1 (SLC22A6), which is critical for maintaining homeostasis of endogenous substances. This makes OAT1 a good transporter to assay for drug interactions with the kidney. Unfortunately, primary cells lose OAT1 expression in culture, and transiently expressed OAT1 has great variations between production lots, which makes data hard to interpret. In our study, we generated HEK 293T/17 cells that stably overexpress the OAT1 gene driven by the human elongation factor-1 alpha (EF1a) promoter. We validated that the overexpressed OAT1 transporter has normal transport activities by using 5-carboxyfluorescein (5-CF) and para-aminohipurate (PAH; data not shown) uptake assays, and that the uptake can be inhibited by the well-known inhibitors probenecid and novobiocin. Overall, our data has shown that this modified cell line is a very useful in vitro tool for testing regulation of OAT1 membrane transporter activity in kidney cells.

MATERIALS AND METHODS

HEK 293T/17 (ATCC[®] <u>CRL-11268</u>[™]) cells were transfected with a plasmid expressing the full-length sequence of human OAT1 (NM_004790) under the control of the EF1α promoter using TransfeX[™] Transfection reagent (ATCC[®] <u>ACS-4005</u>[™]). Transfected cells were grown under antibiotic selection and individual positive clones were identified and verified by RT-PCR, immunofluorescence, and Sanger sequencing. A clone negative for OAT1 or the parental line was used as a control. Cell culture medium consisted of DMEM (ATCC[®] <u>30-2002</u>[™]) supplemented with 10% FBS (ATCC[®] <u>30-2020</u>[™]). OAT1-HEK transfectants were seeded onto poly-L-ornithine coated chamber slides and cultured at 37°C/5% CO₂. After 24 hours, cells were fixed and stained with a rabbit monoclonal antibody against OAT1 (Abcam) and a goat anti-rabbit 488 secondary antibody then visualized using a fluorescent microscope. Sanger sequencing confirmed that the resulting OAT1 HEK 293T/17 (OAT1-HEK; ATCC[®] <u>CRL-11268G-1</u>[™]) line expressed the complete human OAT1 gene with no mutations (data not shown). Copy number was determined via Droplet Digital[™] PCR (Bio-Rad) to be 6 copies per cell (data not shown).

For the 5-CF uptake assay, OAT1-HEK or control cells were seeded at 10^5 cells/well in black walled 96-well plates. After 24h, the cells were washed three times in warm HBSS (ATCC[®] <u>30-2213</u>[™]) and incubated for 10 minutes at 37°C/5% CO₂. Cells were then incubated with 150 μ M 5-CF (Sigma) for 20 min at 37°C/5% CO₂. After incubation, the reaction was terminated by washing the cells three times with cold

HBSS. Cells were then lysed with M-Per Mam*malian* Protein Extraction Reagent (Thermo Fisher Scientific) and read immediately on a fluorescent plate reader at 490ex/530em or visualized by fluorescent microscopy. For the inhibition assay, OAT1-HEK cells were incubated with 5-CF and either probenecid (Sigma) or novobiocin (Sigma) for 10 minutes in 96-well plates, then uptake was measured as described above.

RESULTS/DISCUSSION:

We confirmed the expression of OAT1 in OAT1-HEK by RT-PCR and immunocytochemistry. An approximate 20-fold increase in OAT1 compared to whole kidney lysates was observed. In addition, immunofluorescence revealed that the cells were positive for membrane-localized OAT1 (Figure 1). We then tested the ability of the OAT1-HEK line to uptake a known substrate of OAT1, 5-CF. We observed an uptake ratio of 23:1 relative to the control HEK cells. Moreover, 90% of the OAT1-HEK cells emitted after treatment with the fluorescent 5-CF (Figure 2). The activity of the OAT1-HEK cells was time- and dose-dependent (Figure 3), and stable over 14 passages (Figure 4). Taken together, these data indicate that the OAT1-HEK cell line stably expressed a functional OAT1 protein.

We then tested the ability of two inhibitors of OAT1-mediated activity, probenecid and novobiocin, to block 5-CF transport into the OAT-HEK cells (Figure 5). We observed that both compounds displayed potent inhibitory effects on 5-CF transport (probenecid IC50=15.9 μ M; novobiocin IC50=8.1 μ M), providing further evidence to the usefulness of these cells in toxicological testing.

CONCLUSIONS:

We engineered HEK293T/17 to overexpress the human OAT1 transporter to provide an in vitro model that better mimics the in vivo environment for renal toxicity studies. The OAT1-HEK cell line stably expressed high levels of OAT1 mRNA and membrane-localized protein. We then demonstrated the functionality of the overexpressed OAT by time- and dose-dependent uptake of the fluorogenic substrate 5-CF.

Two known inhibitors of OAT1 activity, probenecid and novobiocin, generated IC50 values similar to those previously reported.^{1,2} Therefore, OAT1-HEK cells and the 5-CF uptake assay serves as a useful alternative to liquid chromatography-mass spectrometry or radioactive assays for sensitive and high-throughput screening of OAT1 modulators.



Figure 1: Confirmation of OAT1 expression in OAT1-HEK by RT-PCR and immunocytochemistry. (A) RT-PCR revealed a ~20-fold increase in human OAT1 relative to whole kidney lysate, N=3. (B) The RT-PCR product was analyzed by gel electrophoresis and found to be of the expected size of the human OAT1 sequence. (C) Immunofluorescent staining with a human anti-OAT1 antibody in cultured OAT1-HEK cells revealed the cells were ~90% positive for membrane localized OAT1.



Figure 2: OAT1 mediated uptake of fluorogenic substrate in OAT1-HEK cells. (A) OAT1-HEK or control cells were incubated with the fluorogenic substrate 5-CF as described in the materials and methods. Mean RFUs for the control line was <400 while the signal from the OAT1-HEK line was >9000, N=3. (B) This indicated an uptake ratio of 23 relative to the control, N=3 (middle). (C) Fluorescent microscopy revealed that ~90% of OAT1-HEK, but not control cells, exhibited uptake and accumulation of 5-CF.



Figure 3: Time- and concentration-dependent uptake of 5-CF in OAT1-HEK cells. OAT1-HEK cells incubated with 5-CF for the indicated times and concentrations demonstrated a linear increase in RFUs ($r2 \ge 0.99$) up to 30 minutes, N=4.







Figure 5: Dose-dependent inhibition of 5-CF uptake by OAT1 inhibitors probenecid and novobiocin in OAT1-HEK. OAT1-HEK cells were incubated with the indicated compounds as described in the materials and methods and tested for 5-CF uptake. The results were graphed as (A) raw signal or as (B) percentage uptake relative to non-treated control cells, N=3.

REFERENCES:

- 1 Takeda M, et al. Characterization of organic anion transport inhibitors using cells stably expressing human organic anion transporters. Eur J Pharmacol 419(2-3): 113-120, 2001. PubMed: 11426832
- Duan P, You G. Novobiocin is a potent inhibitor for human organic anion transporters. Drug Metab Dispos 37(6): 1203-1210, 2009.
 PubMed: 19282394



©2022 American Type Culture Collection. The ATCC trademark and trade name, and any other trademarks listed in this publication are trademarks owned by the American Type Culture Collection unless indicated otherwise. Abcam is a trademark of ABCAM, PLC. Thermo Fisher is a trademark of Thermo Fisher Scientific, Inc. Sigma and Sigma-Aldrich are trademarks of Sigma-Aldrich Co. LLC. Droplet Digital

and Bio-Rad are trademarks of Bio-Rad Laboratories, Inc.

These products are for laboratory use only. Not for human or diagnostic use. ATCC products may not be resold, modified for resale, used to provide commercial services or to manufacture commercial products without prior ATCC written approval.