

Evaluation of Fluoromethyl-2,2-difluoro-1-(trifluoromethyl)vinyl Ether ("compound A") Effects on Urine Protein Excretion in Rats Using Mass Spectrometry

Poster No.
P27-T

Keling Dong¹, Marjorie S Minkoff¹, Matthew Willetts¹, James E Carlson¹, Jeffrey D Miller¹, Evan. D. Kharasch²

¹Applied Biosystems, 500 Old Connecticut Path, Framingham, MA, United States; ²Department of Anesthesiology, Washington University, St. Louis, MO, United States.

ABSTRACT

Fluoromethyl-2,2-difluoro-1-(trifluoromethyl)vinyl ether (FDVE or "compound A"), a haloalkene degradant of the volatile anesthetic sevoflurane, is nephrotoxic in rats. FDVE bioactivation mediates the toxicity, but the molecular and cellular mechanisms of toxicification are unknown. FDVE caused rapid and brisk changes in kidney gene expression, providing potential insights into mechanisms of toxicity, and potential biomarkers for nephrotoxicity[1]. Nevertheless, it is unknown whether gene expression changes are reflected in protein expression, or whether such tissue changes would be reflected in excreted urine proteins. This investigation was to evaluate FDVE effects on urine protein excretion using mass spectrometry and 8-plex iTRAQ® reagents for relative quantitation. Results demonstrate that FDVE causes certain alterations in urine protein/peptide excretion. Multiple components were differentially expressed in a time-dependent manner.

INTRODUCTION

Fluoromethyl-2,2-difluoro-1-(trifluoromethyl)vinyl ether (FDVE or "compound A"), a haloalkene degradant of the volatile anesthetic sevoflurane, is nephrotoxic in rats. FDVE bioactivation mediates the toxicity, but the molecular and cellular mechanisms of toxicification are unknown. FDVE caused rapid and brisk changes in kidney gene expression, providing potential insights into mechanisms of toxicity, and potential biomarkers for nephrotoxicity[1]. Nevertheless, it is unknown whether gene expression changes are reflected in protein expression, or whether such tissue changes would be reflected in excreted urine proteins. This investigation was to evaluate FDVE effects on urine protein excretion using mass spectrometry and 8-plex iTRAQ Reagents for relative quantitation.

MATERIALS AND METHODS

After Animal Use Committee approval, Male Fisher 344 rats (250-300g) housed in individual metabolic cages received a single intraperitoneal injection of 0.25 mmol/kg FDVE, and all urine was collected daily for one week, as described previously[2]. Equal volumes of 6 replicate time points were pooled to create assay time point samples. 150 µL of each pool was brought to 2mL PBS. Each sample was applied to an anti-HSA column (POROS® Affinity Depletion Cartridges). The flow-through (albumin depleted) was desalted on a POROS R150 column. The protein was eluted and dried. Samples were then reconstituted in 1M TEAB. 50µg of each sample (from day 0 to day 7) was processed with the 8-plex iTRAQ® reagents according to manufacturer's instructions. The iTRAQ® reagent labeled sample was then subjected to strong cation exchange chromatography separation and nine fractions were collected. Six of the fractions were analyzed using LC-MALDI on the 4800 MALDI TOF/TOF™ Analyzer (AB/MDS SCIEX). MS and MS/MS data were processed and searched against the IPI rat database (ipi.RAT.v3.26.fasta) using ProteinPilot™ Software (ABSciex).

RESULTS

Figure 1. Ion exchange chromatography fractionation

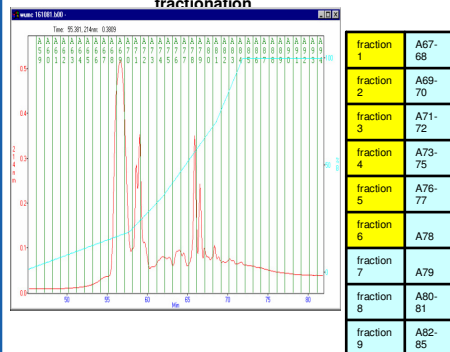
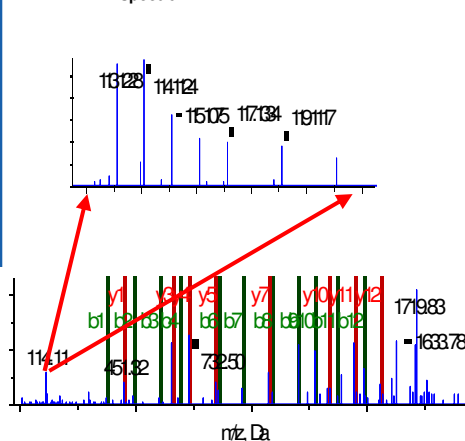


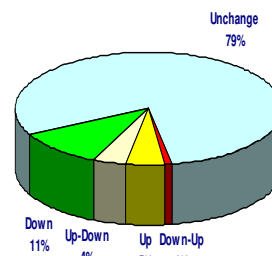
Figure 2. Example of MSMS spectrum.



iTRAQ labeled peptide fragmentation spectrum is shown.

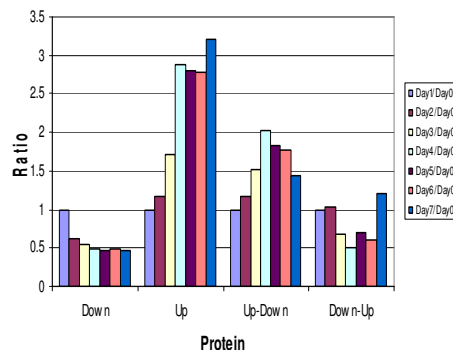
ADLSGITEDAPLK[IT8]
Alpha-1-antiproteinase precursor

Figure 3. Protein expression pie chart



In this experiment there are 231 proteins identified from the FDVE treated rat urine samples. As would be expected, the majority of the identified proteins (79%) show no change in protein expression levels after FDVE treatment. The proteins which do show expression level changes are categorized into 4 groups: There are 11% showing downward trend; 5% upward; 4% up first then down; and 1% down first and then up (see Figure 4 for examples).

Figure 4. Examples of rat urine protein expression level changes after FDVE treatment.



Example	Protein
Down	Haptoglobin precursor
Up	Protein-tyrosine phosphatase LC-PTP
Up-Down	Neuroigin-3 precursor
Down-n-Up	Alpha-2-macroglobulin precursor

There are generally 4 types of trends in the protein level changes: 1) protein level decreases after initial FDVE treatment and slowly levels off; 2) protein level increases after initial FDVE treatment and reaches maximum at the end of sampling; 3) protein level increases initially after FDVE treatment then slowly decreases after reaches maximum; 4) protein level decreases initially after FDVE treatment and slowly levels off then turns upward.

Table 1. List of identified proteins with expression level change after FDVE treatment

Id	Accession #	Name	Trend	U	Accession #	Name	Trend
#6	gpiP1745A1AT	Alpha-1-antiproteinase precursor	Up	#66	gpiP385SC2FU	Cathepsamin precursor	Down
#7	gpiP1234ITRF	Serotransferrin precursor	Up	#67	gpiP1211C2G	Carbonic dehydratase precursor	Down
#8	gpiP2005H4EMO	Hemopexin precursor	Up	#71	gpiQ213N4E9F	Echin-nitin-nitin-binding protein precursor	Down
#12	gpiP2400FETUA	Alpha-2-HS-glycoprotein precursor	Up	#80	gpiP2084E8E2G	Transferrin precursor	Down
#13	gpiP0882ZNTZ	T-lysozyme 2 precursor	Up	#85	gpiQ240M7KFA	Keratin heavy chain isoform 5A	Down
#16	gpiP0906C9P6	Collagenase proteinase inhibitor 6 precursor	Up	#171	gpiQ240M7KFA	Neurokinin B receptor precursor	Down
#17	gpiQ240M7KFA	Fetuin-B precursor	Up	#173	gpiQ240M7KFA	Neurokinin B receptor precursor	Down
#20	gpiP0273TTHY	Transferrin precursor	Up	#184	gpiP2484S7H7	Fibrinogen precursor	Down
#27	gpiP0884HPT	Haptoglobin precursor	Up	#204	gpiP0884HPT	Fibrinogen precursor	Down
#34	gpiP0285SPVIA	Parvalbumin alpha	Up	#204	gpiP0884HPT	Fibrinogen precursor	Down
#35	gpiP0428VTD8	Vitamin D-binding protein precursor	Up	#206	gpiQ20302E2P	EGF precursor 2	Down
#41	gpiP0270TALU	Serum albumin precursor	Up	#211	gpiP0270TALU	Transferrin precursor	Down
#41	gpiP2653AFAM	Albumin precursor	Up	#211	gpiP0270TALU	Transferrin precursor	Down
#43	gpiP0502G2TA1	Glutathione S-transferase alpha-1	Up	#214	gpiP4403RED1	Double-stranded RNA-specific esterase 1	Up-Down
#50	gpiQ2855SPA3M	Serine protease inhibitor AS1 precursor	Up	#240	gpiP288984MP	Serum amyloid P-component precursor	Up-Down
#53	gpiQ2855SPA3M	Serine protease inhibitor AS1 precursor	Up	#240	gpiP288984MP	Serum amyloid P-component precursor	Up-Down
#53	gpiP2764A1AG	Alpha-1-acid glycoprotein precursor	Up	#275	gpiQ2855SPA3M	Transferrin precursor	Up-Down
#59	gpiP2076G2CA	g gamma-2A chain C region	Up	#276	gpiP2076G2CA	Enoyl-CoA hydratase precursor	Up-Down
#73	gpiP288984MP	Cytochrome c, somatic	Up	#276	gpiP2076G2CA	Enoyl-CoA hydratase precursor	Up-Down
#78	gpiP1048MNT1	T-annexin 1 precursor	Up	#281	gpiP2484S7H7	Fibrinogen precursor	Up-Down
#87	gpiP11223THD	Thrombospondin 2 precursor	Up	#281	gpiP2484S7H7	Fibrinogen precursor	Up-Down
#102	gpiQ2484E8E2G	Haptoglobin precursor	Up	#284	gpiP288984MP	Mucin-like associated protein 1B	Up-Down
#120	gpiQ28271EED1	Urokinase-type plasminogen activator	Up	#284	gpiP288984MP	Mucin-like associated protein 1B	Up-Down
#185	gpiQ28271EED1	Zinc finger and homeobox domain 1	Up	#284	gpiP288984MP	Mucin-like associated protein 1B	Up-Down
#185	gpiQ28271EED1	Zinc finger and homeobox domain 1	Up	#284	gpiP288984MP	Mucin-like associated protein 1B	Up-Down
#189	gpiQ28271EED1	Protein Olfact12 homolog	Up	#289	gpiQ28271EED1	Protein Olfact12 homolog	Down-Up

CONCLUSIONS

- The results obtained demonstrate/suggest? that FDVE causes certain alterations in urine protein/peptide excretion.
- Multiple components were differentially expressed in a time-dependent manner. Excretion of several endogenously increased proteins was rapidly decreased by FDVE.
- Other proteins showed increased excretion following FDVE, and then gradually decreased to pre-dose levels.
- Excretion of a third set of proteins, minimally or not detectable in controls, was upregulated following FDVE.
- With the 8-plex iTRAQ® Reagents, relative protein excretion levels can be determined quantitatively, demonstrating that it is an ideal tool for time-course studies.

REFERENCES

- E. D. Kharasch, J. L. Schroeder, T. Bammler, R. Beyer, and S. Srinouprachanh, Toxicological Sciences, 90, 419-431 (2006)
- P. Shaffels, J. L. Schroeder, T. G. Altuntas, H. D. Liggitt, and E. D. Kharasch, Chem. Res. Toxicol. 17, 1177-1189 (2004)

ACKNOWLEDGEMENTS

This project was supported by NIH DK53765

TRADEMARKS/LICENSES

Applied Biosystems is committed to providing the world's leading technology and information for life scientists. Applied Biosystems consists of the Applied Biosystems and Celeris Genomics businesses. Applied Biosystems/MDS SCIEX is a joint venture between Applied Biosystems and MDS Inc. Applied Biosystems, Applied and AB (design) are registered trademarks of Applied Biosystems/MDS SCIEX in the U.S. and/or certain other countries. MIDAS, Paragon, Pro Group, Tempo and ProteinPilot are trademarks and Q TRAP and NanoSpray are registered trademarks of Applied Biosystems/MDS SCIEX. Information subject to change without notice. © 2007 Applied Biosystems and MDS Inc. All rights reserved.