

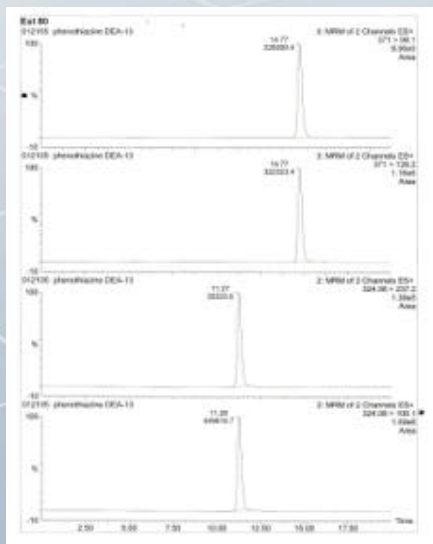
## Drug analysis in animal tissues

### Applications

Within the context of drug analysis in the toxicology field, the responsibility of target molecules in death is investigated. New protocols for lysing matrix are developed to improve the conditions of grinding and the recovery of the molecule regarding constraints of time, cross-contamination and volume of sample.

### Material and method

	Brain	Liver	Heart	Lung	Muscle
<b>Sample</b>	0.5g	0.5g	1g	0.5g	0.3g
<b>Buffer</b>	500µl H <sub>2</sub> O	500µl H <sub>2</sub> O	800µl H <sub>2</sub> O	500µl H <sub>2</sub> O	600µl H <sub>2</sub> O
<b>Kit</b>	ceramic beads 03961CK28	ceramic beads 03961CK28	ceramic beads 03961CK28	ceramic beads 03961CK28	metallic beads 03961MK28
<b>Protocol</b>	Precellys 24 1 cycle 6500 rpm 23s	Precellys 24 1 cycle 6500 rpm 23s	Precellys 24 2 cycles 6500 rpm 23s	Precellys 24 2 cycles 6500 rpm 23s	Precellys 24 2 cycles 6500 rpm 50s
<b>Analysis</b>	LC-MSMS	LC-MSMS	LC-MSMS	LC-MSMS	LC-MSMS



Chromatograms obtained from liver sample spiked with 80 ng/ml of thioridazine and cyamemazine

### Results in collaboration with the Laboratory of Pharmacology - CHU Bordeaux

- time saving compared with manual grinding
- easy to use
- no cross-contamination
- Analysis of liver sample by LC-MSMS :  
Thioridazine recovery 61% and  
cyamemazine recovery 73%