



Figure 2. Calibration curve of the target compounds with concentrations ranging from 0.5 to 80 µg/L.

Conclusion

In this study, an LC/MS/MS method was developed and applied for rapid determination of 44 illicit drugs in urine samples by coupling an LX50 UHPLC with QSight 220 mass spectrometry. The results showed that the method is sensitive and selective with a wide linear range for the studied drugs. Therefore, this method is suitable for fast screening and quantification of these commonly abused drugs in criminal investigations.

References

1. SF/Z JD0107005-2016. Detection of 238 poisons (drugs) in blood and urine.