







Table 4. Method detection limits of TMA using different methods.

Method Name	Sample Volume	Method Detection Limit
Offensive Odor Control Law in Japan	50 L	-
OSHA PV2060	10/20 L	0.2 mg/m <sup>3</sup> (200 µg/m <sup>3</sup> )
HJ 1042-2019	20 L	0.7 µg/m <sup>3</sup>
In This Study	20 L	0.46 µg/m <sup>3</sup>

## Summary

In this study, the determination of TMA in an exhaust pollution source was easily performed utilizing a TurboMatrix HS-40 and Clarus GC/NPD system. The superior detection limit, stability and linearity presented herein ensures a reliable TMA determination.