



ATCC Center for Translational Microbiology

Validation of differentiable molecular barcodes as reference spike-in oligonucleotides for microbiome studies

Tasha M. Santiago-Rodriguez¹, Jessica Schuyler¹, Ron McMillan¹, Gaganpreet Kaur¹, Dev Mittar², John Pace¹, and Raul J. Cano¹ ¹ATCC Center for Translational Microbiology at Institute for Life Science Entrepreneurship, Union, NJ, USA; ²ATCC, Manassas, VA, USA





Figure 5. Taxonomy

Conclusions

- causations.
- significant impact on the results

© 2017 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. Illumina and MiSeq are registered trademarks of Illumina, Inc. PowerSoil and PowerLyzer are registered trademarks of Mo BIO Laboratories, a QIAGEN Company. QIAGEN and QIAamp are registered trademarks of the QIAGEN Group.

phone: 908.737.1922

ATCC-CTM; phone: 908.737.1922; email: info@ilsebio.com







 Results support the value of including reference standards in microbiome analyses to strengthen the validity of findings, correlations, and possible

 Further validation studies are needed to assess the recoverability of the differentiable molecular barcodes and recombinants as in other sample types. • DNA extraction methods, OTU picking strategy, and sequencing depth have a

 The applications of spike-in oligonucleotides can be extended to human and veterinary medicine, personalized medicine, and animal models to study the association of the microbiome in health and disease.