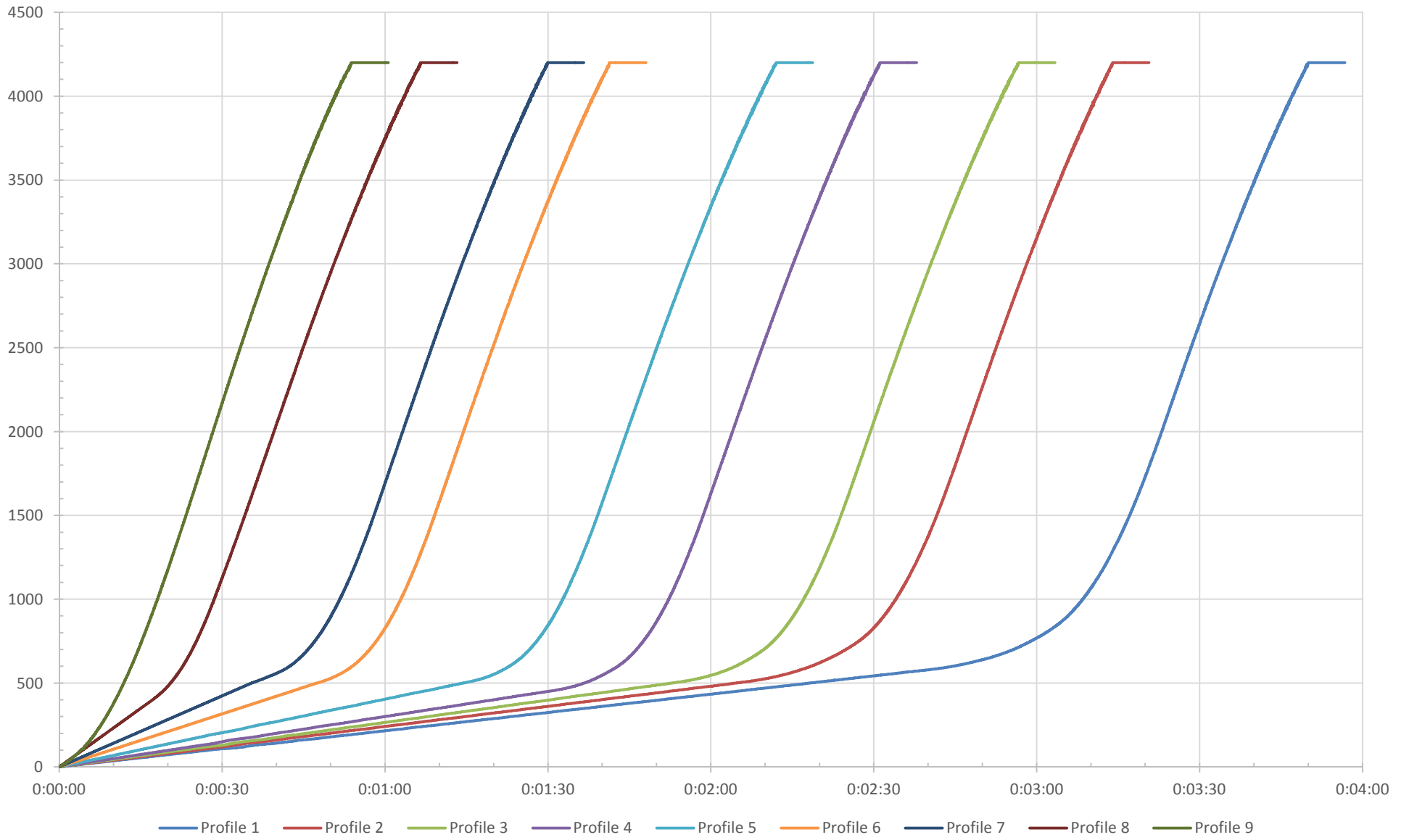


Acceleration profiles TX-1000



Deceleration profiles TX-1000

