



Please provide further information regarding one piece of personal identification according to the following overview:

National IDs / Government IDs	
Country	National IDs / Government IDs
Argentina	DNI Number
Austria	Social Insurance Number (SVNR)
Belgium	ID Card Number
	National Register Number
Brazil	CPF Number
	General Registration (RG) Number
Canada	Social Insurance Number (SIN)
China	Citizen ID Number
Denmark	Central Person Register Number (CPR Nr)
Finland	Personal ID (Henkilötunnus)
France	Social Security ID (INSEE) Number
Germany	Tax ID (Steuer-IdNr)
	Social Insurance Number (SVNR)
Hong Kong	HKID Number
Hungary	ID Card Number
	Social Security (TAJ) Number
India	Aadhaar
Ireland	Personal Public Service Number (PPS no)
Italy	Fiscal Code (CF)
Japan	Individual Number (My Number)
	National Pension Number (Kiso Nenkin Bango)
	Resident (Zairyuu) Card Number
Korea	Resident Registration Number (RRN)



Malaysia	National Identity Number (New IC)
Mexico	Taxpayer (RFC) Number
Netherlands	Personal Number (Burgerservicenr)
Philippines	Social Security System (SSS) Number
Poland	Personal Identity (PESEL) Number
	ID Card Number
Portugal	Social Security Identification Number (NISS)
Puerto Rico	Social Security Number
Singapore	Foreign Identification Number (FIN)
	National Registration Identification Number (NRIC)
Spain	Alien Identification Number (NIE)
	ID Card (DNI) Number
	Social Security Number (NSS)
	Tax Identification Number (NIF)
Sweden	Personal ID Number (Personnummer)
Switzerland	Social Insurance Number (AHV/AVS)
Thailand	ID Card Number
Tunisia	ID Card Number
United Kingdom	National Insurance (NI) Number
United States of America	Social Security Number (SSN)
Vietnam	ID Number