

PUBBLICAZIONE 2022 TRASPARENZA BIO-RAD

I soci devono documentare e rendere pubblici ogni anno attraverso un apposito Modello di Trasparenza che costituisce parte integrante del presente Codice , i trasferimenti di valore effettuati direttamente o indirettamente ai Professionisti del Settore Sanitario, alle Organizzazioni Sanitarie e alle Terze Parti.

	Full Name	HCPs: City of Principal Practice HCOs: City where registered	Country of Principal Practice	Principal Practice Address	Donations	Scholarship	Contribution to costs of Events			Contribution to costs of Training			Costs for the participation in training, educational and promotional events on company's products organised by the Members	Fee for service and consultancy		Transfers of Value Research & Development	Total	
							Sponsorship agreements with HCOs/Third Parties appointed by HCOs to manage an event	Registration Fees	Travel & Accomodation	Sponsorship agreements with HCOs/Third Parties appointed by HCOs to manage an event	Registration Fees	Travel & Accomodation		Fees	Related expenses agreed in the fee for service or consultancy contract, including travel e accommodation relevant to the contract			
HCPs (Healthcare Professionals)	INDIVIDUAL NAMED DISCLOSURE - one line per HCP (I.e. all transfers of value during a year for an individual HCP will be summed up: itemization should be available for the individual Recipient or public authorities' consultation only, as appropriate)																	
					NA	NA	NA*	NA*	NA*							NA	0	
					NA	NA	NA*	NA*	NA*							NA	0	
					NA	NA	NA*	NA*	NA*							NA	0	
					NA	NA	NA*	NA*	NA*							NA	0	
					NA	NA	NA*	NA*	NA*							NA	0	
					NA	NA	NA*	NA*	NA*							NA	0	
					NA	NA	NA*	NA*	NA*							NA	0	
					NA	NA	NA*	NA*	NA*							NA	0	
					NA	NA	NA*	NA*	NA*							NA	0	
					NA	NA	NA*	NA*	NA*							NA	0	
					NA	NA	NA*	NA*	NA*							NA	0	
Healthcare Third Parties	AGGREGATE DISCLOSURE - per HCPs																	
	Aggregate amount attributable to transfers of value to such Recipients				NA	NA	NA*	NA*	NA*				303	1 400	816	NA	2 519	
	Number of Recipients in aggregate disclosure				NA	NA	NA*	NA*	NA*							NA	7	
	% of the number of Recipients included in the aggregate disclosure in the total number of Recipients disclosed				NA	NA	NA*	NA*	NA*							NA	100%	
Healthcare Organizations	INDIVIDUAL NAMED DISCLOSURE - one line per HCOs (I.e. all transfers of value during a year for an individual HCOs will be summed up: itemization should be available for the individual Recipient or public authorities' consultation only, as appropriate)																	
	Biomedica S.r.l.	Milano	Italy	via Libero Temolo, 4	Only HCO	NA	1 220			NA	NA	NA	NA			NA	1 220	
	Centro Italiano Congressi CIC Sud S.r.l.	Bari	Italy	via Escrivà, 28	Only HCO	NA	3 660			NA	NA	NA	NA			NA	3 660	
	Ecomeeting S.r.l.	Catanzaro	Italy	via Gradoni Ospedale Civile, 11	Only HCO	NA	4 270			NA	NA	NA	NA			NA	4 270	
	Koeln Parma Exhibitions S.R.L.	Parma	Italy	viale delle Esposizioni 393a	Only HCO	NA	4 800			NA	NA	NA	NA			NA	4 800	
	S.P.E.M.E - Società per la Promozione dell'Educazione Medica	Catania	Italy	piazza Vincenzo Lanza, 3	Only HCO	NA	4 880			NA	NA	NA	NA			NA	4 880	
	Simtipro S.r.l.	Milano	Italy	via Desiderio, 21	Only HCO	NA	20 130			33 550	NA	NA	NA			NA	53 680	
	Vet International S.r.l.	Milano	Italy	via Carlo Farini, 81	Only HCO	NA	2 702			NA	NA	NA	NA			NA	2 702	
Healthcare Organizations	AGGREGATE DISCLOSURE																	
	Aggregate amount attributable to transfers of value to such Recipients				Only Third Parties		NA	NA	NA	NA	NA	NA	NA	NA	NA			
	Number of Recipients in aggregate disclosure				Only Third Parties		NA	NA	NA	NA	NA	NA	NA	NA	NA			
	% of the number of Recipients included in the aggregate disclosure in the total number of Recipients disclosed				Only Third Parties		NA	NA	NA	NA	NA	NA	NA	NA	NA			

* In the case of direct support for the formation of HCPs that they exercise in a private context, it will be necessary to publish the data in individual or aggregate form, depending on whether the HCP has given consent